# Datasheet

V. 3.1

## Cytoskeleton, Inc.

## MemGlow™ 640: Fluorogenic Membrane Probe

A MEMBRIGHT<sup>™</sup> Family probe

Cat. # MG04-02 (2 nmoles)

he Protein Experts

## Lot #:

Upon arrival store at 4°C (desiccated)

See below for storage after reconstitution

## Background

The MemGlow<sup>™</sup> product line consists of bright & non toxic live cell membrane probes. MemGlow<sup>™</sup> fluorogenic probes exhibit ideal microscopy characteristics including high specificity, low background, and simple application. MemGlow<sup>™</sup> 640 has been validated with multiple microscopy techniques including epifluorescent (widefield), confocal, 2-photon, and TIRF<sup>1</sup>. MemGlow<sup>™</sup> has been confirmed to work in fixed cells, fixed tissue, live cells, and other phospholipid membranes such as extracellular vesicles including exosomes<sup>1-3</sup>.

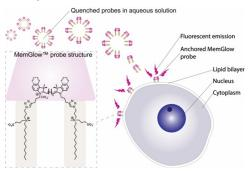


Figure 1. Turn-on mechanism of MemGlow™ probes. MemGlow™ probes are self-quenched nanoparticles until integration with the plasma membrane enables their excitation.

## Material

As measured in DMSO, the absorption max of MemGlow<sup>TM</sup> 640 is 650 nm, with an emission spectra of 673 nm, an extinction coefficient of  $250\times10^3$  cm<sup>-1</sup> M<sup>-1</sup>, and can be visualized using a Cy5<sup>TM</sup> filter set or other suitable filter sets. MemGlow<sup>TM</sup> 640 is supplied as a lyophilized pellet. Avoid contact with MemGlow<sup>TM</sup> by wearing appropriate PPE and dispose of according to local regulations and policies.

#### Storage and Reconstitution

The lyophilized product is stable at 4°C (<10% humidity) for 6 months and should be protected from light. To reconstitute, briefly centrifuge to collect the product at the bottom of the tube. MemGlow<sup>TM</sup> should be reconstituted with 100 µl of anhydrous DMSO to create a 20 µM stock solution for cell imaging or with 10 µl of anhydrous DMSO to create a 200 µM stock suitable for tissue or small organism imaging. After reconstitution the solution should be stored at -20°C where it is stable for 3 months. Once reconstituted, allow product to warm to room temperature before opening tube.

## Online Datasheet Contains MG04-02 (V3.1) and MG04-010 (V3.1)

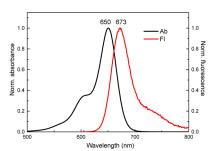


Figure 2. Absorbance and excitation spectra of MemGlow™ 640 diluted in DMSO. Absorbance peak detected at 650 nm and emission peak detected at 673 nm.

## Important Technical Notes

- A. Diluted solutions of MemGlow<sup>™</sup> in aqueous media must be used immediately (<20 sec), as MemGlow<sup>™</sup> will precipitate and/or bind to tube walls.
- B. Serum can reduce MemGlow<sup>™</sup> staining efficiency. When possible MemGlow<sup>™</sup> staining should take place in the absence of serum. Optimally, the imaging cell media is serum-free media, reduced serum media, or PBS. In lieu of serum removal, the concentration of MemGlow<sup>™</sup> should be increased.
- C. Samples incubating in MemGlow<sup>™</sup> solution should be protected from light.
- D. MemGlow™ is non-toxic and live cells can be returned to normal cell media following labeling, and relabeled after 3-4 days.
- E. The localization of MemGlow™ to lipid bilayers is easy to achieve with this product; however, differences in cell morphology and microscope technology, e.g., confocal vs. epifluorescence, will influence the visualization of MemGlow™ (see Figure 3).
- F. When co-labeling with antibodies that require permeabilization limit the concentration of Triton-X to 0.1%.
- G. MemGlow<sup>™</sup> is fully compatible with 4% paraformaldehyde (PFA); however, 4% PFA partially permeabilizes the cell membrane so internalization of probes should be expected.
- H. For tissues and small organisms an initial labeling concentration of 2 μM is recommended. For cell culture an initial labeling concentration of 20-200 nM is recommended depending on application (Table 1).
- Homogeneity of tissue labeling can be optimized with a longer incubation at 4°C rather than relatively brief incubations at room temperature; however, both approaches can label plasma membranes.

cytoskeleton.com

Phone: (303) 322.2254 Fax: (303) 322.2257 Customer Service: cserve@cytoskeleton.com Technical Support: tservice@cytoskeleton.com

## Cytoskeleton, Inc.

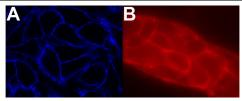


Figure 3. KB (A) or HEK293 (B) cell monolayer labeled with MemGlow™ 640. A) Laser scanning confocal imaging of live KB cells labeled with MemGlow™ 640. Laser excitation was set to 635 nm with emission collected between 640 and 800. B) Widefield fluorescent imaging of live HEK293 labeled with MemGlow™ 640. Cells were imaged as described in **Application 1** methods. Cells were visualized with a Cy5<sup>™</sup> filter set, a digital CCD camera, and 40x objective.

|                                      | Live<br>cells | Fixed cells | Tissue or small<br>organisms |
|--------------------------------------|---------------|-------------|------------------------------|
| Epifluorescent<br>working            | 100           | 100         | 2000                         |
| Confocal<br>working<br>solution (nM) | 20            | 20          | 2000                         |

 Table 1. Recommended initial concentrations. Optimal conditions for efficient labeling should be determined for each cell line and application.

Application 1: Labeling the plasma membrane of live cells in culture.

## Reagents

- 1. MemGlow™ 640 (Cat. # MG04).
- 2. Semi-confluent Tib-71 or HEK293 cells grown in a chamber slide.
- Imaging medias: PBS, serum-free media or reduced serum media.

## Equipment

- Fluorescent microscope with a Cy5<sup>™</sup> excitation filter at 630 +/-20 nm and emission filter at 680 +/-20 nm for MemGlow<sup>™</sup> 640.
- 2. Digital camera.

### Method

- Cells should be seeded onto imaging-appropriate glass or plastics and grown according to cell line requirements to semi-confluency.
- Remove any cell culture media from your cells and replace with the media used for imaging (e.g., serum-free media). Do not allow the cells to dry.
- Prepare the probe solution by diluting 5 µl of 20 µM MemGlow<sup>™</sup> stock in 1 mL imaging media to create a 100 nM working solution or and mix thoroughly. Work quickly (<20 secs) as the probes will begin to aggregate reducing labeling efficiency.
- 4. Add diluted probe solution to cells by replacing the cell media with diluted probe solution until covered. Incubate cells in MemGlow<sup>™</sup> solution for 10 minutes at room temperature. 37°C incubation can be used but will accelerate endocytosis of probes.
- No washing step is required prior to imaging, but can be performed if desired with imaging media.
- Proceed with imaging.

Application 2: Labeling the plasma membrane of fixed cells in culture.

#### Reagents

- 1. MemGlow<sup>™</sup> 640 (Cat. # MG04).
- Semi-confluent Tib71 or HEK293 cells grown on acidwashed coverslips.
- Phosphate-buffered saline (PBS, 20 mM potassium phosphate pH 7.4, 150 mM NaCl).
- Fixative solution (4.0 % paraformaldehyde in PBS).
- Glass microscope slide.
- Coverslip sealing solution (clear nail polish).
- 7. EMS Fluoro-Gel mounting media (Cat. # 17985-10)

## Equipment

- Fluorescent microscope with a Cy5<sup>™</sup> excitation filter at 630 +/-20 nm and emission filter at 680 +/-20 nm for MemGlow<sup>™</sup> 640.
- Digital camera.

## Method

- Cells should be seeded onto imaging-appropriate glass or plastics and grown according to cell line requirements to semi-confluency.
- Remove cell media and wash cells 1X-2X with PBS.
- Fix cells for 10-15 minutes at room temperature with 4% paraformaldehyde (PFA).
- 4. Remove excess PFA by washing cells with PBS 3X.
- (Optional) If co-labeling, permeabilization can be performed at this point. Add 0.1% Triton-X 100 in PBS followed by the primary and secondary antibody protocol according to supplier.
- Prepare the probe solution by diluting 5 µl of 20 µM MemGlow™ stock in 1 ml PBS to create a 100 nM working solution or and mix thoroughly. Work quickly (<20 secs) as the probes will begin to aggregate reducing labeling efficiency.
- Incubate cells in MemGlow<sup>™</sup> solution for 10 minutes at room temperature.
- Remove MemGlow<sup>™</sup> solution and wash cells with PBS 1X-2X.
- 9. If desired place mounting media onto microscope slide.
- 10. Apply cover slip cell-side down onto mounting media or microscope slide.
- 11. If desired apply coverslip sealing solution according to manufacturers directions.
- 12. Proceed with imaging.

## Product Citations/Related Products

- Collot, M. et al. MemBright: A Family of Fluorescent Membrane Probes for Advanced Cellular Imaging and Neuroscience. Cell Chem. Biol. 26, 600-614.e7 (2019).
- Hyenne, V. et al. Studying the Fate of Tumor Extracellular Vesicles at High Spatiotemporal Resolution Using the Zebrafish Embryo. Dev. Cell 48, 554-572.e7 (2019).
- Collot, M., Boutant, E., Lehmann, M. & Klymchenko, A. S. BODIPY with Tuned Amphiphilicity as a Fluorogenic Plasma Membrane Probe. Bioconjug. Chem. 30, 192–199 (2019).

For the latest protocols, citations and related products please visit www.cytoskeleton.com/memglow.

MEMBRIGHT<sup>™</sup> is a trademark of CNRS/UNISTRA of France.

## cytoskeleton.com

# Datasheet

## Cytoskeleton, Inc.

## MemGlow™ 640: Fluorogenic Membrane Probe

A MEMBRIGHT<sup>™</sup> Family probe

```
Cat. # MG04-10 (10 nmoles)
```

he Protein Experts

## Lot #:

Upon arrival store at 4°C (desiccated)

## See below for storage after reconstitution

## Background

The MemGlow<sup>™</sup> product line consists of bright & non toxic live cell membrane probes. MemGlow<sup>™</sup> fluorogenic probes exhibit ideal microscopy characteristics including high specificity, low background, and simple application. MemGlow<sup>™</sup> 640 has been validated with multiple microscopy techniques including epifluorescent (widefield), confocal, 2-photon, and TIRF<sup>1</sup>. MemGlow<sup>™</sup> has been confirmed to work in fixed cells, fixed tissue, live cells, and other phospholipid membranes such as extracellular vesicles including exosmes<sup>1-3</sup>.

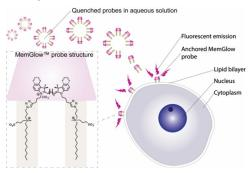


Figure 1. Turn-on mechanism of MemGlow™ probes. MemGlow™ probes are self-quenched nanoparticles until integration with the plasma membrane enables their excitation.

#### Material

As measured in DMSO, the absorption max of MemGlow<sup>TM</sup> 640 is 650 nm, with an emission spectra of 673 nm, an extinction coefficient of  $250\times10^3$  cm<sup>-1</sup> M<sup>-1</sup>, and can be visualized using a Cy5<sup>TM</sup> filter set or other suitable filter sets. MemGlow<sup>TM</sup> 640 is supplied as a lyophilized pellet. Avoid contact with MemGlow<sup>TM</sup> by wearing appropriate PPE and dispose of according to local regulations and policies.

#### Storage and Reconstitution

The lyophilized product is stable at 4°C (<10% humidity) for 6 months and should be protected from light. To reconstitute, briefly centrifuge to collect the product at the bottom of the tube. MemGlow<sup>TM</sup> should be reconstituted with 500 µl of anhydrous DMSO to create a 20 µM stock solution for cell imaging or with 50 µl of anhydrous DMSO to create a 200 µM stock suitable for tissue or small organism imaging. After reconstitution the solution should be stored at -20°C where it is stable for 3 months. Once reconstituted, allow product to warm to room temperature before opening tube.

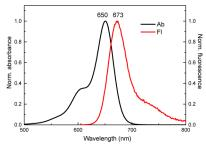


Figure 2. Absorbance and excitation spectra of MemGlow™ 640 diluted in DMSO. Absorbance peak detected at 650 nm and emission peak detected at 673 nm.

## Important Technical Notes

- A. Diluted solutions of MemGlow<sup>™</sup> in aqueous media must be used immediately (<20 sec), as MemGlow<sup>™</sup> will precipitate and/or bind to tube walls.
- B. Serum can reduce MemGlow<sup>™</sup> staining efficiency. When possible MemGlow<sup>™</sup> staining should take place in the absence of serum. Optimally, the imaging cell media is serum-free media, reduced serum media, or PBS. In lieu of serum removal, the concentration of MemGlow<sup>™</sup> should be increased.
- C. Samples incubating in MemGlow<sup>™</sup> solution should be protected from light.
- D. MemGlow™ is non-toxic and live cells can be returned to normal cell media following labeling, and relabeled after 3-4 days.
- E. The localization of MemGlow™ to lipid bilayers is easy to achieve with this product; however, differences in cell morphology and microscope technology, e.g., confocal vs. epifluorescence, will influence the visualization of MemGlow™ (see Figure 3).
- F. When co-labeling with antibodies that require permeabilization limit the concentration of Triton-X to 0.1%.
- G. MemGlow<sup>™</sup> is fully compatible with 4% paraformaldehyde (PFA); however, 4% PFA partially permeabilizes the cell membrane so internalization of probes should be expected.
- H. For tissues and small organisms an initial labeling concentration of 2 μM is recommended. For cell culture an initial labeling concentration of 20-200 nM is recommended depending on application (Table 1).
- Homogeneity of tissue labeling can be optimized with a longer incubation at 4°C rather than relatively brief incubations at room temperature; however, both approaches can label plasma membranes.

cytoskeleton.com

Phone: (303) 322.2254 Fax: (303) 322.2257 Customer Service: cserve@cytoskeleton.com Technical Support: tservice@cytoskeleton.com

## Cytoskeleton, Inc.

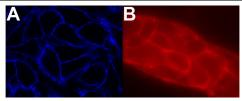


Figure 3. KB (A) or HEK293 (B) cell monolayer labeled with MemGlow™ 640. A) Laser scanning confocal imaging of live KB cells labeled with MemGlow™ 640. Laser excitation was set to 635 nm with emission collected between 640 and 800. B) Widefield fluorescent imaging of live HEK293 labeled with MemGlow™ 640. Cells were imaged as described in **Application 1** methods. Cells were visualized with a Cy5<sup>™</sup> filter set, a digital CCD camera, and 40x objective.

|                                      | Live<br>cells | Fixed cells | Tissue or small<br>organisms |
|--------------------------------------|---------------|-------------|------------------------------|
| Epifluorescent<br>working            | 100           | 100         | 2000                         |
| Confocal<br>working<br>solution (nM) | 20            | 20          | 2000                         |

 Table 1. Recommended initial concentrations. Optimal conditions for efficient labeling should be determined for each cell line and application.

Application 1: Labeling the plasma membrane of live cells in culture.

## Reagents

- 1. MemGlow™ 640 (Cat. # MG04).
- 2. Semi-confluent Tib-71 or HEK293 cells grown in a chamber slide.
- Imaging medias: PBS, serum-free media or reduced serum media.

## Equipment

- Fluorescent microscope with a Cy5<sup>™</sup> excitation filter at 630 +/-20 nm and emission filter at 680 +/-20 nm for MemGlow<sup>™</sup> 640.
- 2. Digital camera.

### Method

- Cells should be seeded onto imaging-appropriate glass or plastics and grown according to cell line requirements to semi-confluency.
- Remove any cell culture media from your cells and replace with the media used for imaging (e.g., serum-free media). Do not allow the cells to dry.
- Prepare the probe solution by diluting 5 µl of 20 µM MemGlow<sup>™</sup> stock in 1 mL imaging media to create a 100 nM working solution or and mix thoroughly. Work quickly (<20 secs) as the probes will begin to aggregate reducing labeling efficiency.
- 4. Add diluted probe solution to cells by replacing the cell media with diluted probe solution until covered. Incubate cells in MemGlow<sup>™</sup> solution for 10 minutes at room temperature. 37°C incubation can be used but will accelerate endocytosis of probes.
- No washing step is required prior to imaging, but can be performed if desired with imaging media.
- Proceed with imaging.

Application 2: Labeling the plasma membrane of fixed cells in culture.

#### Reagents

- 1. MemGlow<sup>™</sup> 640 (Cat. # MG04).
- Semi-confluent Tib71 or HEK293 cells grown on acidwashed coverslips.
- Phosphate-buffered saline (PBS, 20 mM potassium phosphate pH 7.4, 150 mM NaCl).
- Fixative solution (4.0 % paraformaldehyde in PBS).
- Glass microscope slide.
- Coverslip sealing solution (clear nail polish).
- 7. EMS Fluoro-Gel mounting media (Cat. # 17985-10)

## Equipment

- Fluorescent microscope with a Cy5<sup>™</sup> excitation filter at 630 +/-20 nm and emission filter at 680 +/-20 nm for MemGlow<sup>™</sup> 640.
- Digital camera.

## Method

- Cells should be seeded onto imaging-appropriate glass or plastics and grown according to cell line requirements to semi-confluency.
- Remove cell media and wash cells 1X-2X with PBS.
- Fix cells for 10-15 minutes at room temperature with 4% paraformaldehyde (PFA).
- 4. Remove excess PFA by washing cells with PBS 3X.
- (Optional) If co-labeling, permeabilization can be performed at this point. Add 0.1% Triton-X 100 in PBS followed by the primary and secondary antibody protocol according to supplier.
- Prepare the probe solution by diluting 5 µl of 20 µM MemGlow™ stock in 1 ml PBS to create a 100 nM working solution or and mix thoroughly. Work quickly (<20 secs) as the probes will begin to aggregate reducing labeling efficiency.
- Incubate cells in MemGlow<sup>™</sup> solution for 10 minutes at room temperature.
- Remove MemGlow<sup>™</sup> solution and wash cells with PBS 1X-2X.
- 9. If desired place mounting media onto microscope slide.
- 10. Apply cover slip cell-side down onto mounting media or microscope slide.
- 11. If desired apply coverslip sealing solution according to manufacturers directions.
- 12. Proceed with imaging.

## Product Citations/Related Products

- Collot, M. et al. MemBright: A Family of Fluorescent Membrane Probes for Advanced Cellular Imaging and Neuroscience. Cell Chem. Biol. 26, 600-614.e7 (2019).
- Hyenne, V. et al. Studying the Fate of Tumor Extracellular Vesicles at High Spatiotemporal Resolution Using the Zebrafish Embryo. Dev. Cell 48, 554-572.e7 (2019).
- Collot, M., Boutant, E., Lehmann, M. & Klymchenko, A. S. BODIPY with Tuned Amphiphilicity as a Fluorogenic Plasma Membrane Probe. Bioconjug. Chem. 30, 192–199 (2019).

For the latest protocols, citations and related products please visit www.cytoskeleton.com/memglow.

MEMBRIGHT<sup>™</sup> is a trademark of CNRS/UNISTRA of France.

## cytoskeleton.com