

Cat #	MG15-10
Name	eFM488 Enhanced fluorogenic membrane probe
Amount	10 nmoles
Lot #	051

Test	Specification	Lot results
Appearance	Orange film	Pass
Absorption max. in DMSO	483 nm	Pass
Cell membrane staining test (1 μ M eFM488)	Membrane staining of live A431 cells by fluorescence microscopy using FITC filter set (excitation 450-480 nm and emission 515-555). No nuclear staining	Pass
Concentration	When resuspended in 100 μ l of DMSO the concentration is 100-125 μ M	124 μ M
QC release date	2/9/26	
Approved by	K. Middleton	