

Cat #	MG16-10
Name	eFM560 Enhanced fluorogenic membrane probe
Amount	10 nmoles
Lot #	001

Test	Specification	Lot results
Appearance	Purple film	Pass
Absorption max. in DMSO	557 nm	Pass
Cell membrane staining test (1 $\mu$ M eFM560)	Membrane staining of live A431 cells by fluorescence microscopy using filter set (excitation 562 nm and emission 624). No nuclear staining	Pass
Concentration	When resuspended in 100 $\mu$ l of DMSO the concentration is 100-125 $\mu$ M	118 $\mu$ M
QC release date	2/9/26	
Approved by	K. Middleton	