

Cat #	TL488M
Name	Tubulin protein (HiLyte 488 labeled) from Porcine Brain
Amount	20 µg
Lot #	037

Test	Specification	Lot results
Appearance	Light orange lyophilized powder	Pass
Purity by Coomassie Gel	≥99%	99%
Protein Quantitation with Precision Red Protein Assay Reagent (Cytoskeleton Cat. # ADV02)	20 – 30 µg per tube	27.3 µg
Biological Activity Assay: Polymerization Assay	>75% Polymerization	80%
Labeling Stoichiometry: HiLyte 488 dyes per tubulin protein molecule	0.5 – 1.0 dyes per protein molecule	0.7
QC release date	6/26/26	
Approved by	K. Middleton	