

Cat #	TL670M
Name	Tubulin protein (HiLyte 647 labeled) from Porcine Brain
Amount	20 µg
Lot #	041

Test	Specification	Lot results
Appearance	Light blue 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 µg
Biological Activity Assay: Polymerization Assay	>75% Polymerization	78%
Labeling Stoichiometry: HiLyte 647 dyes per tubulin protein molecule	0.1 – 0.3 dyes per protein molecule	0.12
QC release date	9/29/25	



Kim Middleton
 QC Manager