

MICAL1 protein (Human)
 RedoxCH p45 domain, 6xHis
 Store at 4°C (desiccated)

Catalog # MIC01
 MIC01-XL

Product Description: Human MICAL1 protein (MIC01) [accession # Q8TDZ2 (MICA1_HUMAN)] has been produced and purified from a bacterial expression system. The recombinant protein is N-terminally 6x Histidine tagged and is comprised of the Redox and CH domains (aa: 1-612) of the full length protein. MIC01 has an approximate molecular weight of 60 kDa. MIC01 is supplied as a yellow lyophilized powder.

Amount: MIC01 50 µg
 MIC01-XL 1 mg

Lot: 001

TEST	SPECIFICATION		LOT RESULTS
Appearance	Yellow lyophilized powder		Yellow lyophilized powder
Purity by Coomassie Gel	>90%		>90%
Protein Quantitation with Precision Red Protein Assay	MIC01	50-60µg	53.3 µg
Reagent (Cytoskeleton Cat. # ADV02)	MIC01-XL	1.0-1.2 mg	na
Biological Activity Assay #1: Subtilisin cleavage	<10% cleavage of oxidized Actin. >90% cleavage of native actin		<10% >90%
Biological Activity Assay #2: Spin-down polymerization assay	40-50% polymer at 0.1 mg/ml 70-80% polymer at 0.4 mg/ml		40% 80%
QC Release Date	April 17 th 2019		



Kim Middleton
 QC Manager