

Pyrene Labeled Actin Protein (Mical-oxidized)

Rabbit skeletal

Store at 4°C (desiccated)

Catalog # MXAP95
MXAP95-XL

Product Description: Pyrene labeled rabbit skeletal muscle actin protein (MICAL-oxidized) (Cat. # MXAP95) has been enzymatically oxidized at methionines 44 and 47 (b-actin nomenclature) with the MICAL flavoprotein monooxygenase protein (Cat. # MIC01). MICAL-oxidized (pyrene labeled) rabbit muscle actin has an approximate molecular weight of 43 kDa and is supplied as a white lyophilized powder.

Amount: MXAP95 250 µg
MXAP95-XL 1 mg

Lot: 001

TEST	SPECIFICATION	LOT RESULTS
Appearance	White lyophilized powder	White lyophilized powder
Purity by Coomassie Gel	>95%	>95%
Protein Quantitation with Precision Red Protein Assay Reagent (Cytoskeleton Cat. # ADV02)	MXAP95 250-275µg MXAP95-XL 1.0-1.2 mg	261 µg 1.04 mg
Biological Activity Assay #1: Subtilisin cleavage	<10% cleavage of Mical1 oxidized actin. >75% cleavage of native actin	<10% >75%
Biological Activity Assay #2: Pyrene actin polymerization assay (fluorescence: ex 360 nm, em 405 nm) of oxidized pyrene actin compared to native pyrene actin	At 0.05 mg/ml actin: oxidized pyrene actin will not polymerize (<10% above baseline) At 0.1 mg/ml actin: oxidized actin will polymerize at 40-50% the efficiency of native pyrene actin	<10% 50%
QC Release Date	May 29 th 2019	



Kim Middleton



Certificate of Analysis



QC Manager

To Place Order Tel: (303) 322-2254, Fax: (303) 322-2257, E-mail: cserve@cytoskeleton.com
For Technical Assistance E-mail: tservice@cytoskeleton.com. Cytoskeleton, Inc. Denver, CO. USA
WWW.CYTOSKELETON.COM