

ProQinase™ MRCL3

myosin light chain 12A

Recombinant Protein Kinase Substrate

HGNC Symbol: MYL12A

Synonyms: MLC-2B, MLCB, MRCL3, MYL2B, RLC

Product No.: 1575-0000-1

Lot: 001

Description: MRCL3, full length, amino acids M₁-D₁₇₁ (as in [NCBI/Protein](#) entry NP_006462.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in E.coli

Theoretical MW_{Fusion Protein}: 48,697 Da

Expression host: E.coli

Purification: GST-Affinity Chromatography

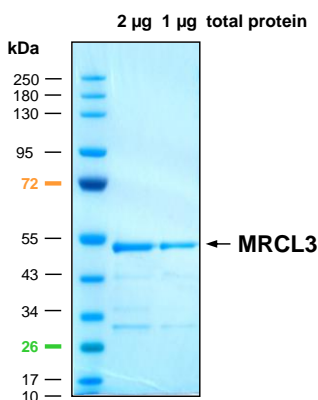
Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 10 % glycerol

Storage temperature: -80°C

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

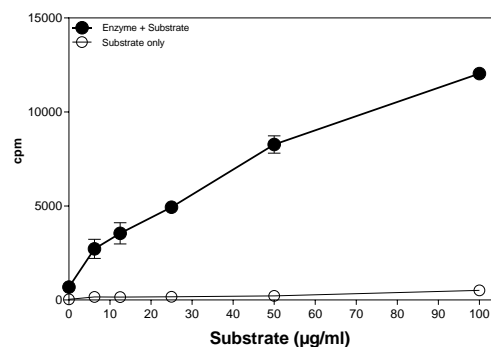
Protein concentration: 0.854 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

MRCL3 Lot 001: Coomassie stain



Phosphorylation of MRCL3 by STK17B

Radiometric filter binding assay



Assay conditions:

70 mM HEPES-NaOH, pH 7.5
3 mM MgCl₂
3 mM MnCl₂
3 µM Na-orthovanadate
1.2 mM DTT
50 µg/ml PEG_{20,000}
ATP: 1 µM
Substrate: variable concentration
Kinase: 2 µg/ml
MSFC membrane (Millipore)

Sequence information

GST-MRCL3 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGQATF	GGGDHPPKSD	PMG HHHHHG	RDS LEVLFGQ	240
241	PLAMGARGRM	SSKRTKTKTK	KRPQRATSNV	FAMFDQSQIQ	EFKEAFNMID	QNRDGFIDKE	300
301	DLHDMLASLG	KNPTDEYLDA	MMNEAPGPIN	FTMFLTMFGE	KLNGTDPEDV	IRNAFACFDE	360
361	EATGTIQEDY	LRELLTTMGD	RFTDEEVDEL	YREAPIDKKG	NFNIEFTRI	LKHGAKDKDD	420

1-218: GST **Red:** HIS6-tag **Green:** 3C cleavage site **blue:** MRCL3

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in human or animals.
 © European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH