

ProQinase™ CFL1

cofilin 1

Recombinant Protein Kinase Substrate

HGNC Symbol: CFL1

Synonyms: CFL

Product No.: 1753-0000-1

Lot: 001

Description: CFL1, full length, amino acids M₁-L₁₆₆ (as in [NCBI/Protein](#) entry NP_005498.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in E.coli

Theoretical MW_{Fusion Protein}: 47,007 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

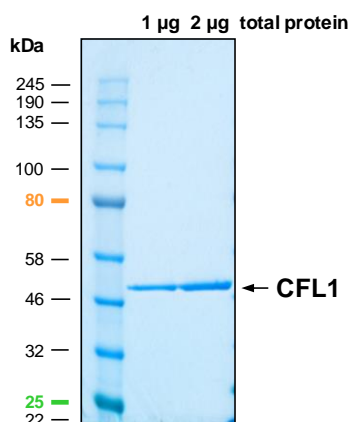
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 2.2 µg/µl

(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

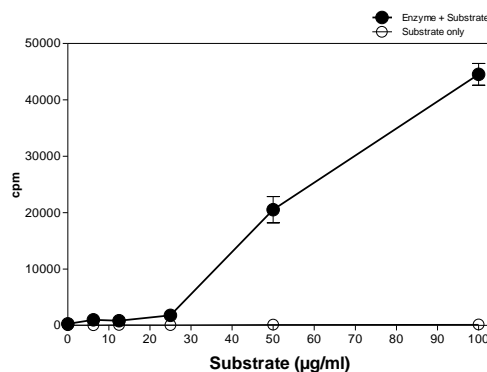
CFL1 Lot 001:

Coomassie stain



Phosphorylation of CFL1 by LIMK1

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: 1 µg/ml
MSFC membrane (Millipore)

Field of application:

CFL1 has been validated for use in radiometric in-vitro kinase activity assays. It was not validated for use in ATP-consumption based kinase activity assays.

ProQinase™ CFL1

Product No.: 1753-0000-1

GST-CFL1 Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGQ	240
241	PLAMVMASGV	AVSDGVIKVF	NDMKVRKSST	PEEVKKRKA	VLFLSEDKK	NIILEEGKEI	300
301	LVGDVGQTV	DPYATFVKML	PDKDCRYALY	DATYETKESK	KEDLVFIFWA	PESAPLKSKM	360
361	IYASSKDAIK	KKLTGIKHEL	QANCYEEVKD	RCTLAEKLGG	SAVISLEGKP	L	420

1-218: GST Red: HIS6-tag Green: 3C cleavage site blue: CFL1

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