

ProQinase™ CDCA8

cell division cycle associated 8

Recombinant Protein Kinase Substrate

HGNC Symbol: CDCA8

Synonyms: borealin, FLJ12042, MESRGP, BOR, DasraB

Product No.: 1773-0000-1

Lot: 001

Description: CDCA8, full length, amino acids M₁-K₂₈₀ (as in [NCBI/Protein](#) entry NP_060571.1), N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in E.coli

Theoretical MW_{Fusion Protein}: 60,226 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

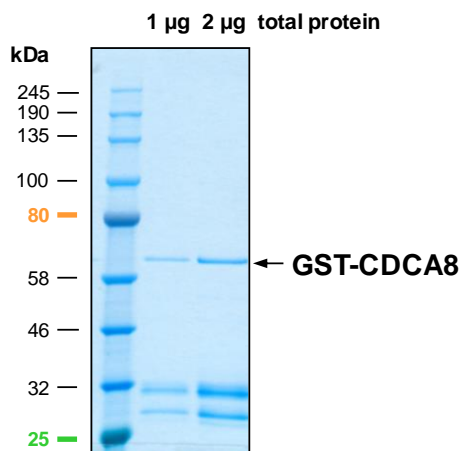
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 1.332 µg/µl

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

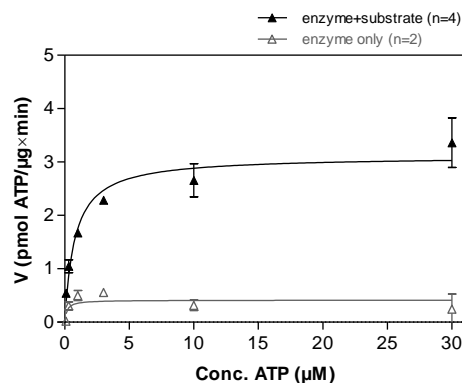
Substrate Lot 001:

Coomassie stain



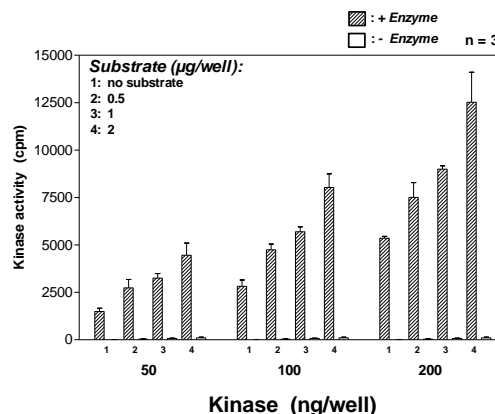
Phosphorylation of CDCA8 by TTK

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: CDCA8 20 µg/ml
 Kinase TTK: 2 µg/ml
 MSFC membrane (Millipore)

³³PanQinase Assay (Flashplate based)



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}
 1 µM ATP
 1 % (v/v) DMSO

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

ProQinase™ CDCA8

Product No.: 1773-0000-1

CDCA8 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFKDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQG WQATF	GGGDHPPKSD	PMGHHHHHG	RDSLEVLFGG	240
241	PLAMGARGRM	APRKGSSRVA	KTNSLRRRKL	ASFLKDFDRE	VEIRIKQIES	DRQNLLKEVD	300
301	NLYNIEILRL	PKALREMNWL	DYFALGGNKQ	ALEEAATADL	DITEINKLTA	EAIQTPLKSA	360
361	KTRKVIQVDE	MIVEEEEEEE	NERKNLQTAR	VKRCPPSKKR	TQSIQKGKGG	KRSSRANTVT	420
421	PAVGRLEVSM	VKPTPGLTPR	FDSRVFKTPG	LRTPAAGERI	YNISGN GSPL	ADSKEIFLTV	480
481	PVGGGESLRL	LASDLQRHSI	AQLDPEALGN	IKKLSNRLAQ	ICSSIRTHK		540

1-218: GST **Green**: 3C cleavage site **blue**: CDCA8

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