

ProQinase™ RBER-CDC25tide

Generic Recombinant Protein Kinase Substrate

HGNC Symbol: n/a

Synonyms: n/a

Product No.: 0590-0000-1

Lot: 012

Description: Artificial fusion protein consisting of an N-terminal GST-tag separated by a Thrombin cleavage site from a fragment of the human retinoblastoma protein RB1, amino acids S₇₇₃-K₉₂₈ (as in [NCBI/Protein](#) entry NP_000312.2) followed by 11 Arg residues (ER) and a peptide sequence (ISDELMDATFADQEAQ; CDC25tide). Expressed in E.coli.

Theoretical MW_{Fusion Protein}: 47,608 Da

Expression host: E.coli

Purification: GST-Affinity Chromatography

Field of application:

JNK1 K55R/K56R has been validated for use in radiometric in-vitro kinase activity assays.

It is not recommended for use in ATP-consumption based kinase activity assays.

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

Storage temperature: -80°C

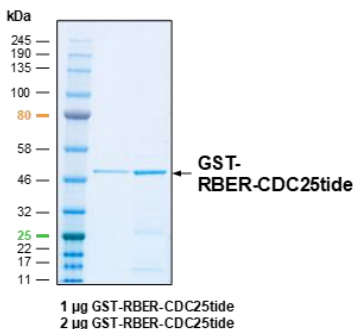
For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10 µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 1.274 µg/µl

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

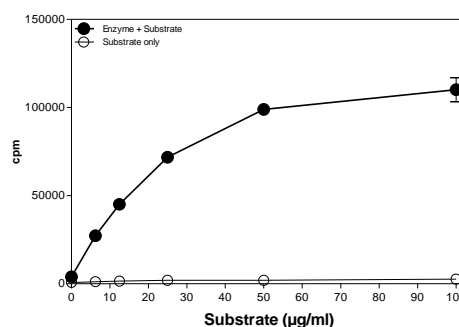
RBER-CDC25tide Lot 012:

Coomassie stain



Phosphorylation of RBER-CDC25tide by PLK1

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)

ProQinase™ RBER-CDC25tide

Product No.: 0590-0000-1

RBER-CDC25tide Recombinant Fusion Protein Amino Acid Sequence

1	MSPILGYWKI	KGLVQPTRLL	LEYLEEKYEE	HLIERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLPEML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAIPQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	LVPRGSPPEFS	TRPPTLSPIP	240
241	HIPRSPYKFP	SSPLRIPGGN	IYISPLKSPY	KISEGLPTPT	KMTPRSRIIV	SIGESFGTSE	300
301	KFQKINQMC	NSDRVLKRS	EGSNPPKPLK	KLRFDIEGSD	EADGSKHLP	ESKFQQLAE	360
361	MTSTRTRMQ	QKMNDSDTS	NKEEKRRRR	RRRRRISDE	LMDATFADQ	EAK	420

1-218: GST **Pink**: Thrombin cleavage site **blue**: RB1 fragment **Green**: R₁₁-sequence **boxed**: CDC25tide sequence