

ProQinase™ RBER-NTRK3tide

Recombinant Protein Kinase Substrate

Product No.: 1132-0000-1

Lot: 011

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 (VYSTDYRFLNPS), derived from the human NTRK3 protein ([NCBI/Protein](#) entry NP_001012338.1, aa V704-S716).

Theoretical MW_{Fusion Protein}: 47,450 Da

Expression host: E.coli

Purification: GST-Affinity and ion exchange chromatography

ATPase activity: In an ADP-Glo™ assay (Promega) with 10 μM ATP or 30 μM ATP, the ATP → ADP conversion within 30 min is < 1% at a concentration of 100 μg/ml substrate.

Detailed ATPase assay conditions on request

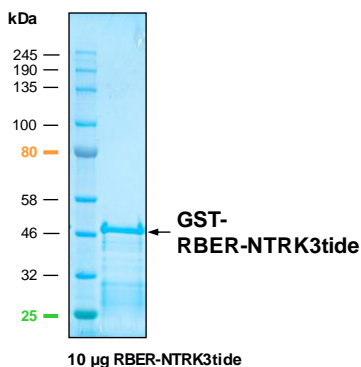
Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20 % glycerol

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!

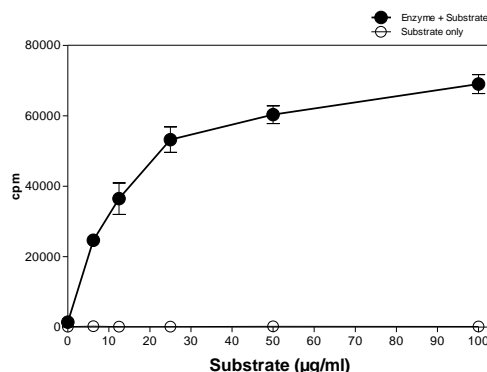
Protein concentration: 2.325 μg/μl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

RBER-NTRK3tide Lot 011:
Coomassie stain



Phosphorylation of RBER-NTRK3tide by TRK-C

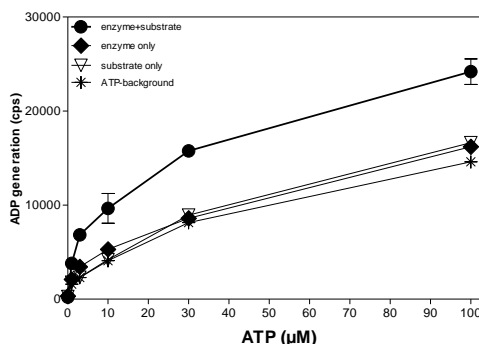
Radiometric filter binding assay



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: 1 μg/ml
MSFC membrane (Millipore)

ADP-Glo™ assay (Promega)



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: variable concentration
1 % (v/v) DMSO
Substrate: 100 μg/ml
Kinase: 1 μg/ml

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™ RBER-NTRK3tide

Product No.: 1132-0000-1

RBER-NTRK3tide 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	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGDHPPKSD	LVPRGSP EFS	TRPPTLSPI P	240
241	HIPRSPYKFP	SSPLRIPGGN	IYISPLKSPY	KISEGLPTPT	KMTPRSRILV	SIGESFGTSE	300
301	KFQKINQ MVC	NSDRVLKRSA	EGSNPPKPLK	KLRFDIEGSD	EADGSKHLP G	ESKFQQLAE	360
361	MTSTRTRMQ K	QKMND SMDTS	NKEEKRRRRR	RRRRR R	VYST	DYRLFNPS	420

1-218: GST **Pink**: Thrombin cleavage site **Green**: R₁₁-sequence **blue**: RB1 fragment **boxed**: NTRK3tide sequence

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