

RBER-NTRK3tide

Recombinant Protein Kinase Substrate

Product No.: 1132-0000-1

Lot: 002

Description: Artificial fusion protein consisting of a 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 (VYSTDYRLFNPS), derived from the human NTRK3 protein (GenBank entry Q 16288, aa V₇₀₄-S₇₁₆). Expressed in E.coli.

Theoretical MW: 47,450 Da

Expression: E. coli

Purification: Affinity chromatography using GSH-agarose, followed by ion exchange chromatography

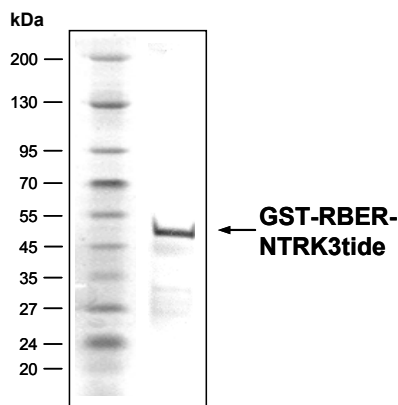
ATPase activity: In an ADP-Glo™ assay (Promega) with 10 µM ATP or 30 µM ATP, the ATP → ADP conversion within 30 min is below 1% at a concentration of 100 µg/ml RBER-NTRK3tide*
*detailed ATPase assay conditions on request

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 1 mM DTT, 15 mM reduced glutathione, 10% glycerol

Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

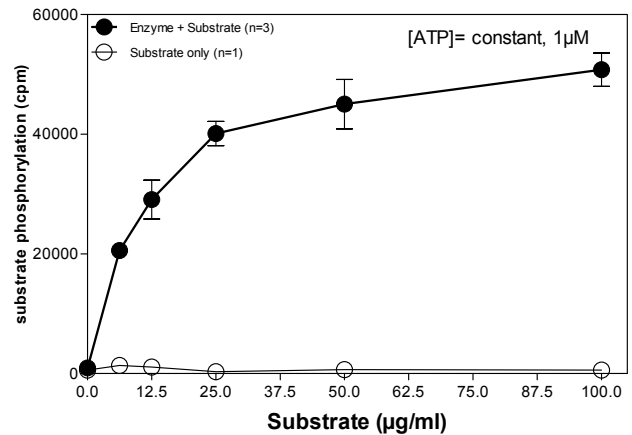
Protein concentration: 1.010 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Coomassie stain:



3.0 µg GST-RBER-NTRK3tide

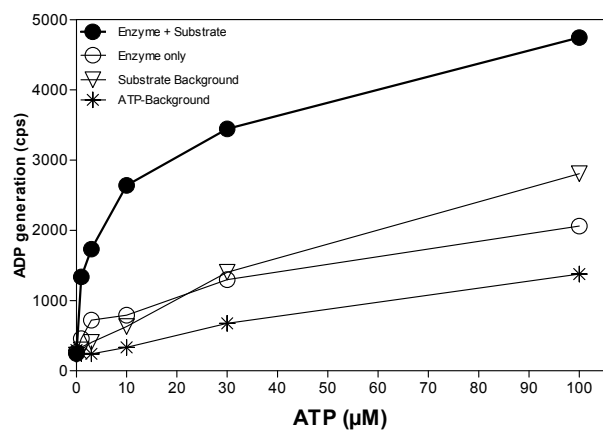
Phosphorylation of RBER-NTRK3tide by the kinase TRK-C (Radiometric filter binding assay):



Assay mixture:

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 RBER-NTRK3tide: variable concentration
TRK-C: 800 ng/ml
MSFC membrane (Millipore)

Phosphorylation of RBER-NTRK3tide by the kinase TRK-C (ADP-Glo™ assay / Promega):



Assay mixture:

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 RBER-NTRK3tide: 100 µg/ml
TRK-C: 800 ng/ml

This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.