

ProQinase™ TESTPRODUKT LOT002

XXX standard example full name

Recombinant Human Active Protein Kinase

HGNC Symbol: KIN1

Synonyms: Kanase, Konase

Product No.: 0000-0000-1

Lot: 000

Description: Human Kinase, C/N/internal fragment, amino acids M₁-X₁₀₀₀ (as in [NCBI/Protein](#) entry NP_000000.0), N-terminal GST-HIS₆ fusion protein with a Thrombin/3C cleavage site, expressed in Sf9 insect cells/E.coli

Product identity: Kinase Lot 000, was confirmed as Kinase by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein:} 00,000 Da

Expression host: Sf9 insect cells/E.coli

Purification: GST-Affinity Chromatography
Immobilized Metal Affinity Chromatography

Activation: This kinase was not activated by special procedures / With KinaseY / in vitro auto activation

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 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 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.000 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

Specific kinase activity (P_i transfer): 0 pmol/µg*min

ATP-K_M: 0 µM

Additional assay technology:

Kinase Lot 000 was also successfully tested by Reaction Biology for the use with the ADP-Glo Kinase assay from Promega
ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

Kinase Lot 000:
Coomassie stain

Kinase Lot 000:
Determination of V_{max} and K_M value for ATP

- Assay conditions:
 - 60 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)
 - Substrate: Example 0 µg/ml
 - Kinase: 0 µg/ml
- Filter binding assay
MSFC/PH membrane (Millipore)

Recombinant Proteins

Sequence information

GST-Kinase Recombinant Fusion Protein Amino Acid Sequence		
1		60
61		120
121		180
181		240
241		300
301		360
361		420
421		480
481		540
541		600
601		660

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **Green**: 3C/TEV cleavage site **blue**: Kinase fragment **boxed**: variation from RefSeq

Kinase wt ¹ Amino Acid Sequence		
1		60
61		120
121		180
181		240
241		300
301		360
361		420
421		480
481		540
541		600
601		660
661		720
721		780
781		840
841		900
901		960
961		1020
1021		1080
1081		1140
1141		1200
1201		1260
1261		1320
1321		1380
1381		1440
1441		1500
1501		1560
1561		1620

blue: kinase sequence expressed in recombinant protein **Red**: variant in recombinant protein

¹[NCBI/Protein](#) accession number NP_000000.0
A100B: SNP variation see [NCBI/dbSNP](#) ID: rs0000000

GST-Kinase Recombinant Fusion Protein Amino Acid Sequence		
1		60
61		120
121		180
181		240
241		300
301		360
361		420
421		480
481		540
541		600
601		660

1-218: GST **Red**: HIS6-tag **Pink**: Thrombin cleavage site **Green**: 3C/TEV cleavage site **blue**: Kinase fragment **boxed**: variation from RefSeq

Kinase wt ² Amino Acid Sequence		
1		60
61		120
121		180
181		240
241		300
301		360
361		420
421		480
481		540
541		600
601		660
661		720
721		780
781		840
841		900
901		960
961		1020
1021		1080
1081		1140
1141		1200
1201		1260
1261		1320
1321		1380
1381		1440
1441		1500
1501		1560
1561		1620

blue: kinase sequence expressed in recombinant protein **Red**: variant in recombinant protein

²[NCBI/Protein](#) accession number NP_000000.0
A100B: SNP variation see [NCBI/dbSNP](#) ID: rs0000000
[HGNC](#) identifier #COFACTOR#: YYYY