

p38-alpha K53A (inactivated)
(mitogen-activated protein kinase 14)

Recombinant Protein Kinase Substrate

Synonyms: MAPK14, CSBP1

Product No.: 0769-0000-1

Lot: 007

Description: Human p38-alpha, full length, amino acids M₁-S₃₆₀ (as in GenBank entry NM_001315)*, mutationally inactivated: K53A, N-terminal HIS₆ fusion protein with a TEV cleavage site, expressed in E.coli
*Sequence may contain documented polymorphisms
Detailed aa-sequence on request

Theoretical MW_{Fusion Protein}: 43,615 Da

Expression: E.coli

Purification: Immobilized Metal Affinity Chromatography followed by a size exclusion chromatography

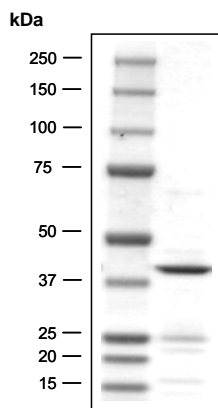
ATPase activity: In an ADP-Glo™ assay (Promega) with 10 µM ATP or 30 µM ATP, the ATP → ADP conversion within 30 min is approx. 1% at a concentration of 100 µg/ml p38-alpha K53A*
*detailed ATPase assay conditions on request

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

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

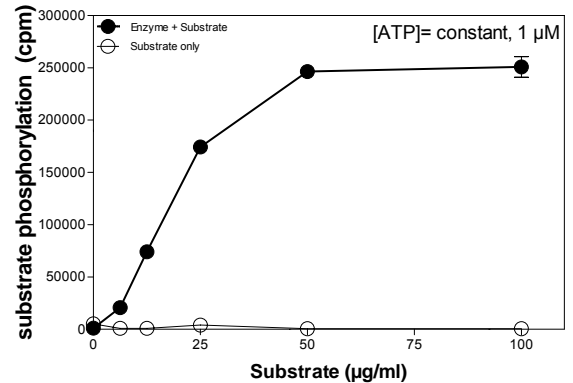
Protein concentration: 0.660 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Coomassie stain:



2 µg p38-alpha K53A

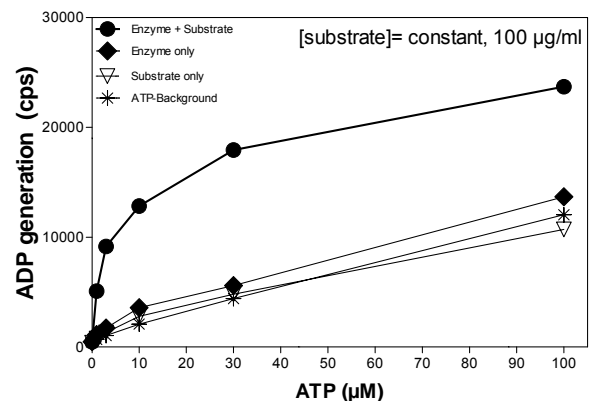
Phosphorylation of p38-alpha K53A by the kinase MKK6 S207D/T211D (Radiometric filter binding assay):



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 (p38-alpha K53A): variable concentration
- MKK6 S207D/T211D: 2.0 µg/ml
- MSFC membrane (Millipore)

Phosphorylation of p38-alpha K53A by the kinase MKK6 S207D/T211D (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 (p38-alpha K53A): 100 µg/ml
- MKK6 S207D/T211D: 2.0 µg/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.