

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

Recombinant Protein Kinase Substrate

Synonyms: MAPK14, CSBP1

Product No.: 0769-0000-1

Lot: 013

**Description:** Human p38-alpha, full length, amino acids M<sub>1</sub>-S<sub>360</sub> (as in NCBI/Protein entry NP\_001306.1), mutationally inactivated: K53A, N-terminal HIS<sub>6</sub> fusion protein with a TEV cleavage site, expressed in E.coli

**Theoretical MW<sub>Fusion Protein</sub>:** 43,615 Da

**Expression:** E.coli

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

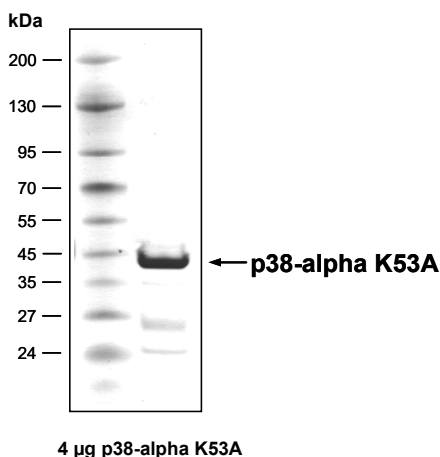
**ATPase activity:** In an ADP-Glo™ assay (Promega) with 10 μM ATP or 30 μM ATP, the ATP → ADP conversion within 30 min is approx. 1% at a concentration of 100 μg/ml p38-alpha K53A\*  
\*detailed ATPase assay conditions on request

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

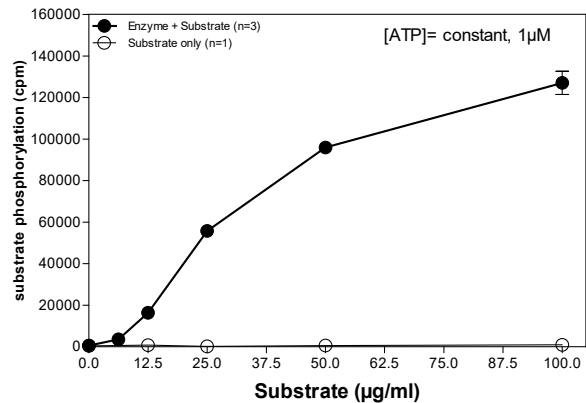
**Storage temperature:** -80°C  
Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.771 μg/μl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Coomassie stain:**



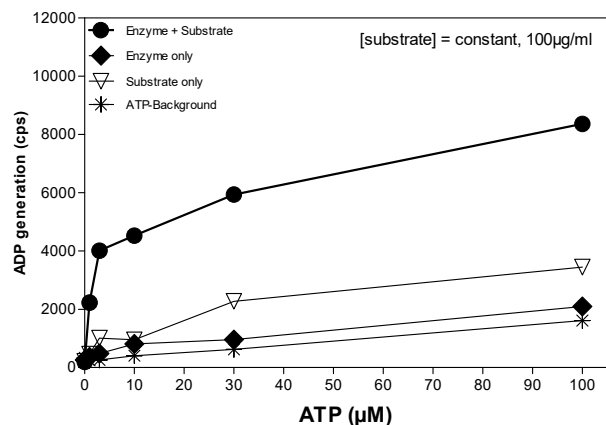
### Phosphorylation of p38-alpha K53A by the kinase MKK6 S207D/T211D (Radiometric filter binding assay):



Assay mixture:

70 mM HEPES-NaOH, pH 7.5  
3 mM MgCl<sub>2</sub>  
3 mM MnCl<sub>2</sub>  
3 μM Na-orthovanadate  
1.2 mM DTT  
50 μg/ml PEG<sub>20,000</sub>  
ATP: 1 μM  
Substrate (p38-alpha K53A): variable concentration  
MKK6 S207D/T211D: 800 ng/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 MgCl<sub>2</sub>  
3 mM MnCl<sub>2</sub>  
3 μM Na-orthovanadate  
1.2 mM DTT  
50 μg/ml PEG<sub>20,000</sub>  
ATP: variable concentration  
1% (v/v) DMSO  
Substrate (p38-alpha K53A): 100 μg/ml  
MKK6 S207D/T211D: 800 ng/ml

Recombinant Proteins