

ProQinase™ TAOK3

TAO kinase 3

Recombinant Human Active Protein Kinase

HGNC Symbol: TAOK3

Synonyms: DPK, JIK, KDS, MAP3K18, hKFC-A

Product No.: 0908-0000-1

Lot: 005

Description: Human TAOK3, full length, amino acids M₁-R₃₉₈ (as in [NCBI/Protein](#) entry NP_057365.2), untagged, expressed in Sf9 insect cells

Product identity: TAOK3 Lot 005, has been verified by mass spectrometry LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 106015 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

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 10µl are not advisable. Avoid repeated freeze-thaw cycles!

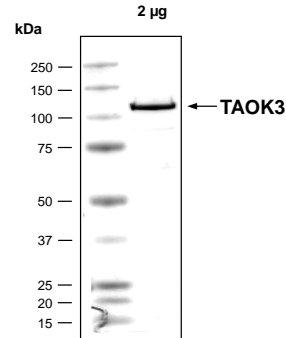
Protein concentration: 0.124 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Biochemical Parameters:

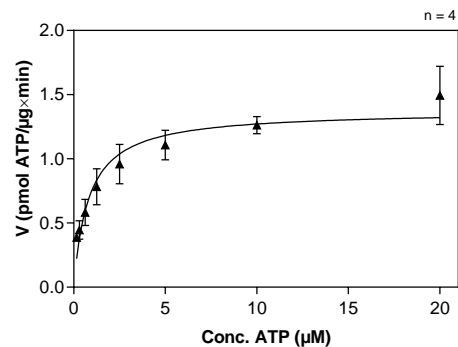
Specific kinase activity (P_i transfer): 1.4 pmol/µg*min

ATP-K_M: 0.81 µM

**TAOK3 Lot 005:
Coomassie stain**



**TAOK3 Lot 005:
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: PKC-derived peptide 2.5 µg/ml
- Kinase: 2 µg/ml

Assay technology:

- Radiometric filter binding assay
- MSPH membrane (96 well plate, Millipore)

Recombinant Proteins

