

## ProQinase™ TAOK3 (TAO kinase 3)

### Recombinant Human Active Protein Kinase

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

**Product No.:** 0908-0000-1

**Lot:** 005

**Description:** Human TAOK3, full length, amino acids M<sub>1</sub>-R<sub>898</sub> (as in GenBank entry NM\_016281.2)\*, untagged, expressed in Sf9 insect cells

\*Sequence may contain documented polymorphisms Detailed aa-sequence on request

**Product identity:** TAOK3 Lot 005 was confirmed as TAOK3 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Protein</sub>:** 106,015 Da

**Expression:** Baculovirus infected Sf9 cells

**Purification:** GST-Affinity Chromatography

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 20% glycerol

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

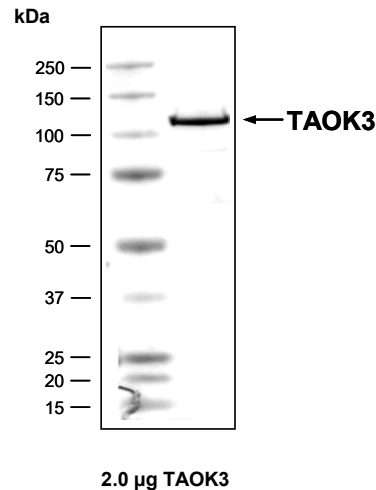
**Protein concentration:** 0.124 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### Determination of K<sub>m</sub> value & Specific activity:

- Assay conditions:
  - 60 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)
  - Substrate: PKC-derived peptide (LRRLSLGLRRLSLGLRRLSLG-RFARKGSLRQKNV), 2.5 µg / ml
  - TAOK3: 2.0 µg / ml
- Filter binding assay
  - MSPH membrane (Millipore)

**Specific activity:** 1 pmol/µg×min

### Coomassie stain:



### Determination of K<sub>m</sub> value for ATP:

