

# FAK aa411-686

focal adhesion kinase

**Synonyms:** PTK2, FADK

**Product No.:** 0480-0000-6

**Lot:** 009

**Description:** Human FAK, internal fragment, amino acids S<sub>411</sub>-Q<sub>686</sub> (as in GenBank entry L13616)\*, untagged, expressed in Sf9 insect cells

\*Sequence may contain documented polymorphisms  
Detailed sequence on request

**Product identity:** FAK, Lot 009, was confirmed as FAK by mass spectroscopy LC-ESI-MS/MS (Protagen AG, Germany)

**Theoretical MW:** 31,807 Da

**Expression:** Baculovirus infected Sf9 cells

**Purification:** One-step affinity purification using Ni-MAC, followed by cleavage of the His<sub>6</sub> tag

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol

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

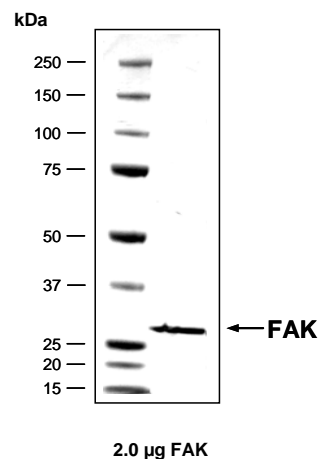
**Protein concentration:** 1.075 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

## Method for 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
  - 2.5 µg / 50 µl PEG<sub>20.000</sub>
  - ATP (variable)
  - Substrate: Poly(Glu:Tyr)<sub>4:1</sub> (Sigma 20K5903), 0.5 µg / 50 µl
  - FAK tr: 200 ng / 50 µl
- Filter binding assay
  - MSFC membrane (Millipore)

**Specific activity:** 18 pmol/µg×min

## Coomassie stain:



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

