

ProQinase™ FAK aa411-686

focal adhesion kinase

Synonyms: PTK2, FADK

Product No.: 0480-0000-6

Lot: 009

Description: Human FAK, internal fragment, amino acids S₄₁₁-Q₆₈₆ (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 His₆ 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!

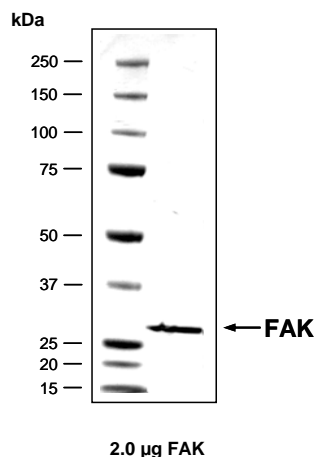
Protein concentration: 1.075 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Method for determination of K_m value & Specific activity:

- Assay conditions:
60 mM HEPES-NaOH, pH 7.5
3 mM MgCl₂
3 mM MnCl₂
3 µM Na-orthovanadate
1.2 mM DTT
2.5 µg / 50 µl PEG_{20,000}
ATP (variable)
Substrate: Poly(Glu:Tyr)_{4:1}
(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_m value for ATP:

