

## ProQinase™ FGF-R1 V561M (fibroblast growth factor receptor 1)

Recombinant Human Active Protein Kinase

Synonyms: b-FGF-R, FLT2

Product No.: 0890-0000-1

Lot: 003

**Description:** Human FGF-R1 V561M, internal fragment, amino acids M<sub>456</sub>-E<sub>765</sub> (as in GenBank entry NM\_000604.2)\*, V561M gatekeeper mutant, untagged, expressed in Sf9 insect cells

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

**Product identity:** FGF-R1 V561M Lot 003 was confirmed as FGF-R1 by mass spectroscopy LC-ESI-MS/MS

**Theoretical MW<sub>Protein</sub>:** 35,899 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!

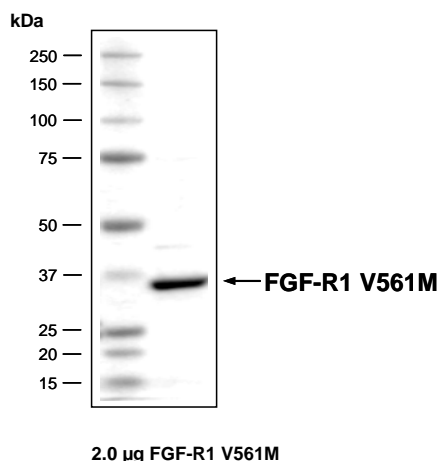
**Protein concentration:** 0.140 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

### 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: Poly(Ala,Glu,Lys,Tyr)<sub>6:2:5:1</sub> (Sigma P-1152), 20 µg / ml
  - FGF-R1 V561M: 0.6 µg / ml
- Filter binding assay  
MSFC membrane (Millipore)

**Specific activity:** 75 pmol/µg×min

### Coomassie stain:



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

