

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 M₄₅₆-E₇₆₅ (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_{Protein}: 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!

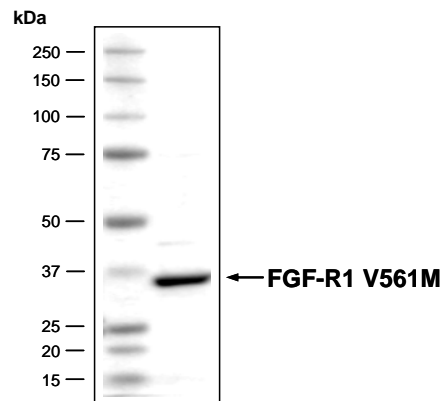
Protein concentration: 0.140 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Determination of K_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
 - 50 µg / ml PEG_{20,000}
 - ATP (variable)
 - Substrate: Poly(Ala,Glu,Lys,Tyr)_{6:2:5:1} (Sigma P-1152), 20 µg / ml
 - FGF-R1 V561M: 0.6 µg / ml
- Filter binding assay
MSFC membrane (Millipore)

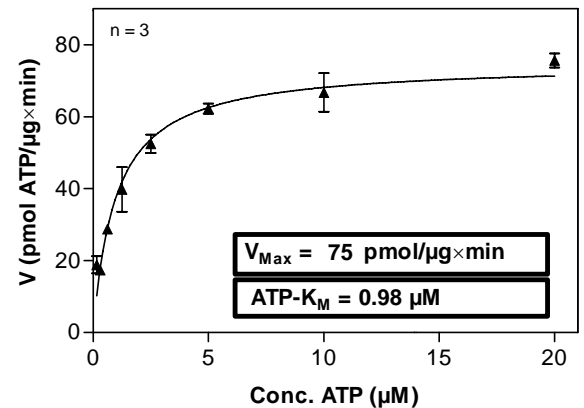
Specific activity: 75 pmol/µg×min

Coomassie stain:



2.0 µg FGF-R1 V561M

Determination of K_m value for ATP:



This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.