

ABL1 T315I

(v-abl Abelson murine leukemia viral oncogene homolog 1)

Recombinant Human Active Protein Kinase

Synonyms: JTK7, p150

Product No.: 0700-0000-1

Lot: 002

Description: Human ABL1, internal fragment, amino acids P₁₁₆-S₅₃₅ (as in GenBank entry NM_005157.3)*, T315I gatekeeper mutant, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

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

Product identity: ABL1 T315I Lot 002, was confirmed as ABL1 by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 78,620 Da

Expression: Baculovirus infected Sf9 cells

Purification: GST-Affinity Chromatography

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol

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

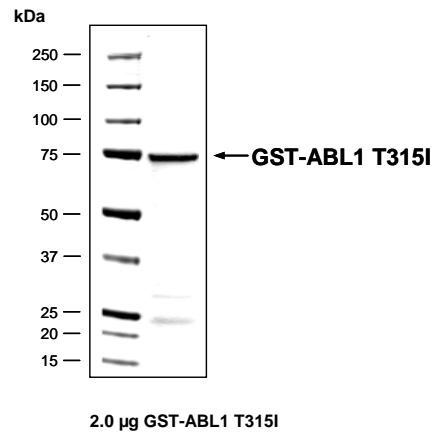
Protein concentration: 0.111 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Determination of K_m value & Specific activity:

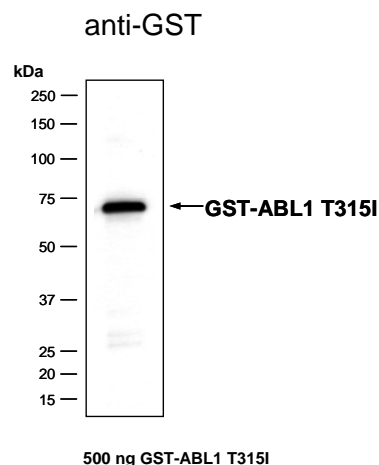
- Assay conditions:
 - 60 mM HEPES-NaOH, pH 7.5
 - 3 mM MgCl₂
 - 3 mM MnCl₂
 - 3 µM Na-orthovanadate
 - 1.2 mM DTT
 - 50 µg / ml PEG_{20,000}
 - ATP (variable)
 - Substrate: Poly(Ala,Glu,Lys,Tyr)_{6:2:5:1} (Sigma P-1152), 20 µg / ml
 - ABL1 T315I: 2.0 µg / ml
- Filter binding assay
MSFC membrane (Millipore)

Specific activity: 36 pmol/µg×min

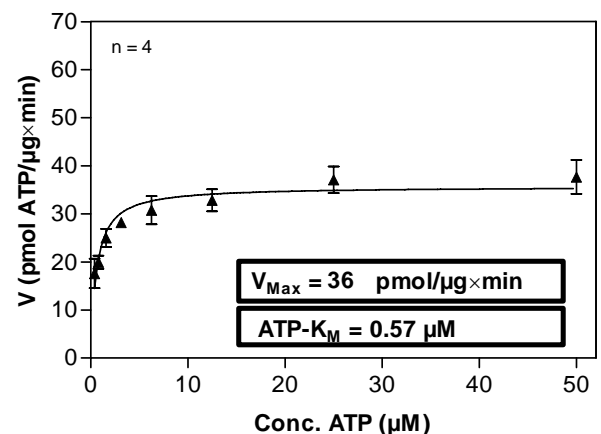
Coomassie stain:



Western blot analysis:



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.