

Data Sheet

Fucokinase

Human, recombinant, N-terminal FLAG-tag
 Catalog # 71077
 Lot#: 130129 Conc.: 0.34 mg/ml

Formulated in: 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween, 3 mM DTT, 20% glycerol, and 80 ng/μL FLAG peptide.

Stability: >6 months at -80 °C

Reference(s):

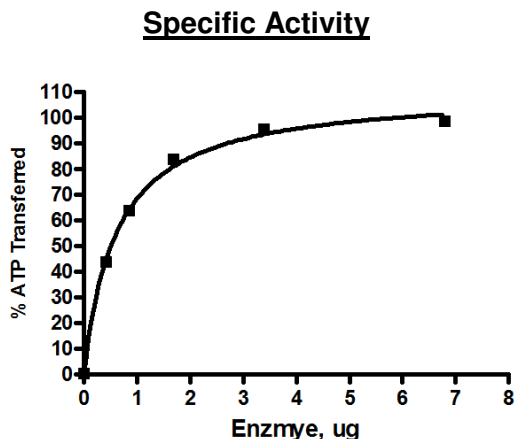
1. Hinderlich, S., *et al.*, *Biochem Biophys Res Commun.* 2002 Jun 14; **294(3)**:650-4.
2. Zeitler, R., *et al.* *J. Enz. Inhib.* 1997 Feb; **11(4)**:265-73.

Description: Human fucokinase (GenBank Accession No. NM_145059), amino-acids 2-1084(end) with N-terminal FLAG-tag, MW = 118.6 kDa, expressed in baculovirus using an Sf9 insect cell expression system.

Specific Activity: 85.3 pmole/min/μg
Assay Conditions: The reaction mixture contained 50 mM Tris-HCl, pH 8.0, 5 mM MgCl₂, 2 mM L-Fucose, 100 μM ATP, and various amount of fucokinase in a final volume of 50 μL. The reaction was incubated at 37 °C for 1 hr, and the amount of ATP consumed was calculated using Kinase-Glo Plus[®] reagent (Promega).

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

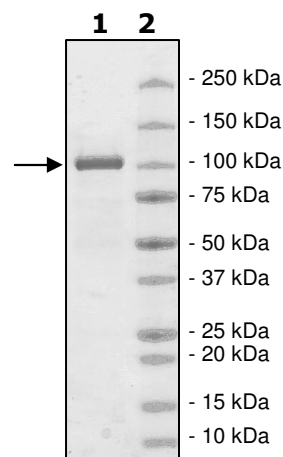
Quality Assurance



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg Fucokinase
Lane 2:
Protein Marker

MW: 118.6 kDa
Purity: ≥95%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.259.5734**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

**MDYKDDDDKEQPKGVDWTVIILTCQYKDSVQVFQRELEVRQKREQIPAGTLLLAVEDPEK
RVGSGGATLNALLVAAEHLSARAGFTVVTSVDLHSAWILILHMGRDFPFDDCGRAFTCLP
VENPEAPVEALVCNLDCLLDIMTYRLGPGSPPGVWVCSTDMLLSVPANPGISWDSFRGA
RVIALPGSPAYAQNHGVYLTDPQGLVLDIYYQGTEAEIQRCVRPDGRVPLVSGVFFSVET
AERLLATHVSPPLDACTYLGLDSGARPVQLSLFFDILHCMAENVTRDFLVGRPPELGQG
DADVAGYLSARAQLWRELRDQPLTMAYVSSGSYSYMTSSASEFLLSLTLPGAPGAQIV
HSQVEEQQLLAAGSSVVSCLLEGPVQLGPGSVLQHCHLQGPIHIGAGCLVTGLDTAHSK
ALHGRELRLDLVQGHHTRLHGSPGHAFTLVGRLD SWERQGAGTYLNVPWSEFFKRTGV
RAWDLWDPETLPAEYCLPSARLFPVLHPSRELGPQDLLWMLDHQEDGGEALRAWRAS
WRLSWEQLQPCLDRAATLASRRDLFFRQALHKARHVLEARQDLSLRPLIWA AVREGCP
GPLLATLDQVAAGAGDPGVAARALACVADVLGCM AEGRGGLRSGPAANPEWMRPF SY
LECGDLAAGVEALAQERDKWLSRPALLVRAARHYEGAGQILIRQAVMSAQHFVSTEQVE
LPGPGQWVVAECPARVDFSGGWS DTPPLAYELGGAVLGLAVRVDGRRPIGARARRIPEP
ELWLAVGPRQDEMTVKIVCRCLADLRDYCQPHAPGALLKAAFICAGIVHVHSELQLSEQL
LRTFGGGFELHTWSELPHGSG LGTSSILAGTALAALQRAAGR VVGTEALIHAVLHLEQVL
TTGGGWQDQVGG LMPGIKVGRSRAQLPLKVEVEEVTVPEGFVQKLN DHLLL VYTGKTRL
ARNLLQDVLRSWYARLP AVVQNAHSLVRQTEECAEGFRQGS LPLL GQCLTSYWEQKKL
MAPGCEPLTVRRMMDVLAPHVHGQSLAGAGGGGFLYLLTKEPQQKEALEAVLAKTEGL
GNYSIHLVEVD TQGLSLKLLGTEASTCCPFP**

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.259.5734**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com