

Data Sheet

FYN A

Human, recombinant, N-terminal Flag tag

Catalog #: 40433

Lot #: 120712 Conc. 0.38 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 80 ng/μl Flag
 peptide, 20% glycerol and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Ban, K. *et al.*, *Blood*, **111** (5), 2409-8 (2008)
2. Hoe, HS. *et al.*, *J. Biol. Chem* **283** (10), 6288-99 (2008)

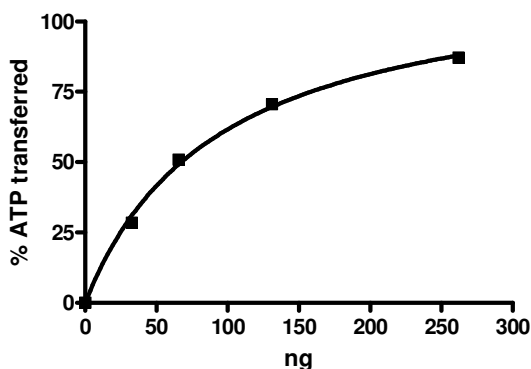
Description: Human FYN A (GenBank Accession No. NM_002037), full length with N-terminal Flag-tag, a.a. 2-537(end), MW= 61 kDa, expressed in Sf9 insect cells.

Specific Activity: 207 pmole/min/μg.
Assay Conditions: Enzyme reaction is conducted in Kinase Buffer (50 mM Tris pH 7.4, 20 mM MgCl₂, 0.1 mg/ml BSA and 1 mM DTT) containing 20 μM ATP and 0.2 mg/ml poly-(Glu-Tyr) for 30 min at 30°C. Enzyme activity is measured using Kinase-Glo Plus™ reagents (Promega).

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

Quality Assurance

Specific Activity

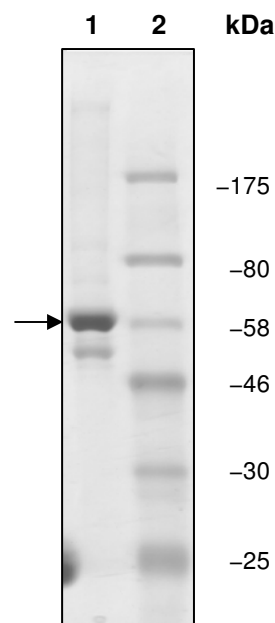


**10% SDS-PAGE
 Coomassie staining**

Lane 1:
 3 μg FYN A

Lane 2:
 Protein Marker

MW: 61kDa
Purity: ≥66%



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 **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



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

Sequence:

MDYKDDDDKGCVQCKDKEATKLT EERDGS LNQSSGYRYGTDPTPQHYP SFGVTSIP
NYNNFHAAGGQGLTVFGGVNSSSHTGTLRTRGGTGVTLFVALYDYEARTEDDLSFHK
GEKFQILNSSEG DWWEARSLTTGETGYIPSNYVAPVDSIQAE EWYFGKLGRKDAERQ
LLSFGNPRGTFLIRESETTKGAYSL SIRDWDDMKGDHVKHYKIRKLDNGGYYITTRAQ
FETLQQLVQHYSERAAAGLCCRLVVPCHKGMPRLTDLSVKT KDVWEIPRESLQLIKRL
GNGQFGEVWMGTWNGNTKVAIKTLKPGTMSPE SFL EEAQIMKCLKHDKLVQLYAVV
SEPIYIVTEYMNKGSLDLDFLKDGEGRALKLPNLVDM AAQVAAGMAYIERMNYIHRDL
RSANILVGNGLICKIADFLARLIEDNEYTARQGAKFPIKWT APEAALYGRFTIKSDVW
SFGILLTELVTKGRVPYPGMNNREVLEQVERGYRMP CPQDCPISLHELMHCWKKDP
EERPTFEYLQSFLEDYFTATEPQYQGENL

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 **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com