



6042 Cornerstone Ct. West, Ste. B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

IRS1-tide

Catalog #: 79519 Lot #: 190731
Size: 0.5 mg @ 1 mg/ml (10X)

Formulation: Proprietary buffer for peptide reconstitution

Stability: At least 6 months at -20°C .

References:

1. Varkondi E. *et al.*, *J of Receptors and Signal Transduction*, 2005; **25(1)**; 45-56.

Description:

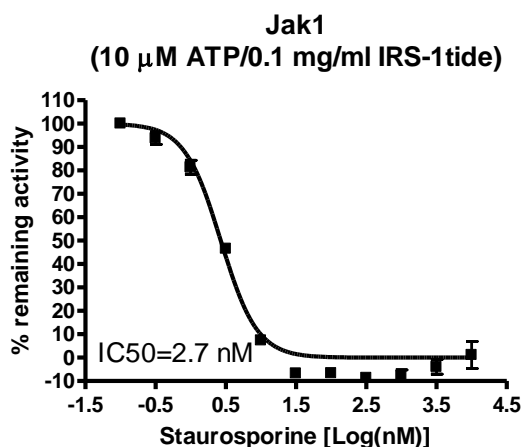
Ready to use Kinase substrate (10X)
IRS1-tide is a peptide derived from the sequence motif of the insulin receptor substrate-1.

Application:

Useful as a substrate for protein tyrosine kinase inhibitor screening including Jak1 Ron kinase, and Tyrosine Kinase 2.

Quality Assurance

IRS-1 tide



Enzyme reaction was performed using Jak1 (Janus Kinase 1) Assay Kit (BPS #79518)

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

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com