

Data Sheet

STAT5 Peptide

Catalog #: 79864 **Lot #:** 191101
Size: 0.5 mg @ 10 mg/ml (50X)

Formulation: Reconstitute in H₂O.

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

References:

1. Saharinen P. *et al.*, *Mol. Cell. Biol.*, 2000;
20(10); 3389-3395.

Description:

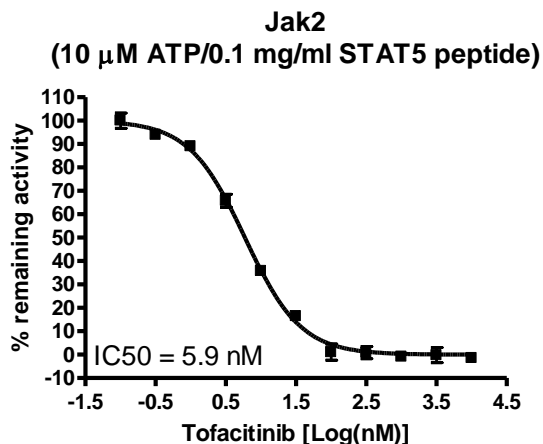
Ready to use Kinase substrate (50X)
STAT5 peptide is a peptide derived from
the sequence motif of the STAT5.

Application:

Useful as a substrate for JAK2 inhibitor
screening.

Quality Assurance

STAT5 Peptide



Enzyme reaction was performed by
following Jak2 (Janus Kinase 2) Assay Kit
(BPS Bioscience #79520) using STAT5
peptide as a substrate.

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