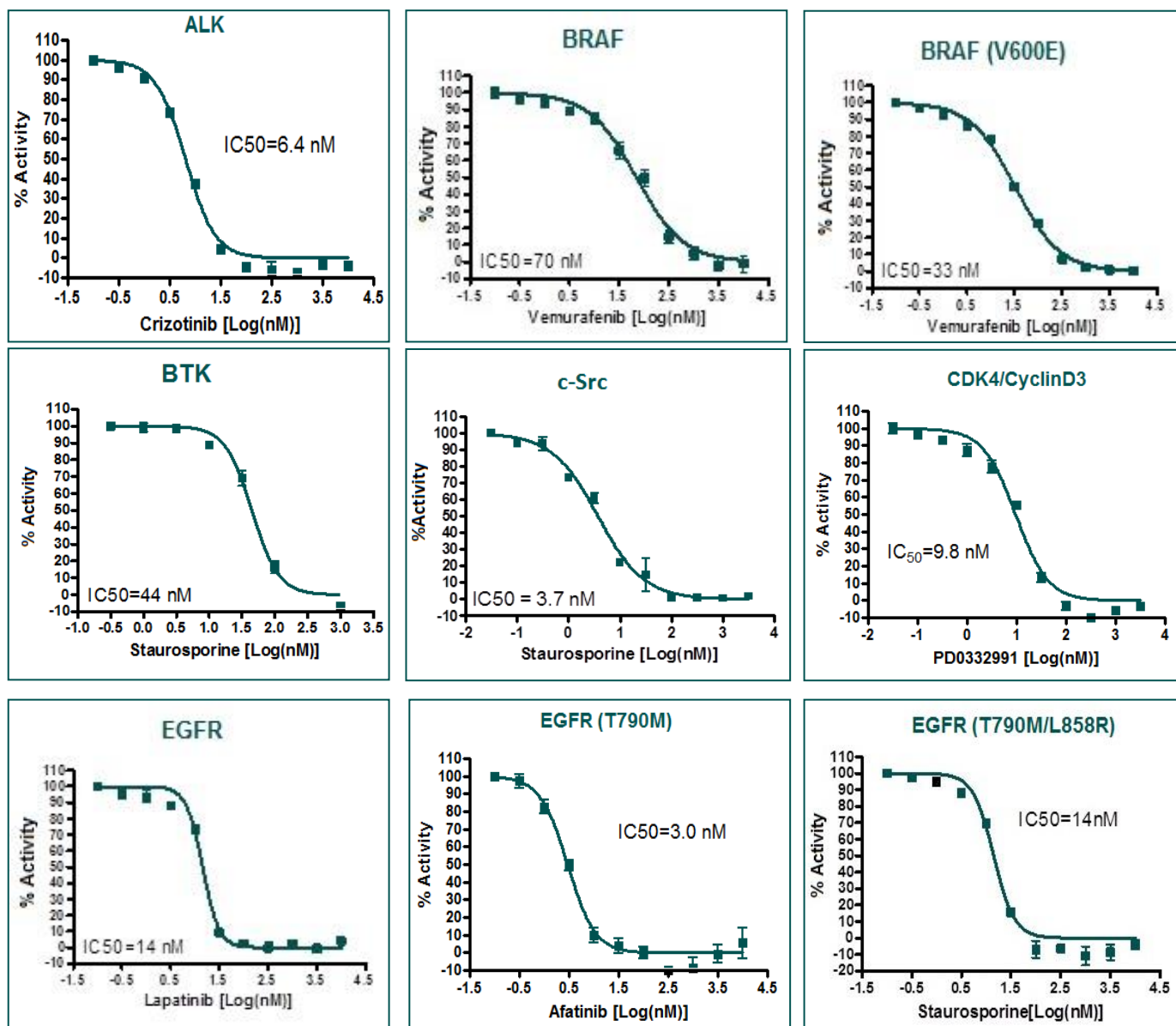


Kinase Screening / Profiling

Determine the IC₅₀ of your compound, or screen against our panel of kinases.
 Contact us for a quote today!

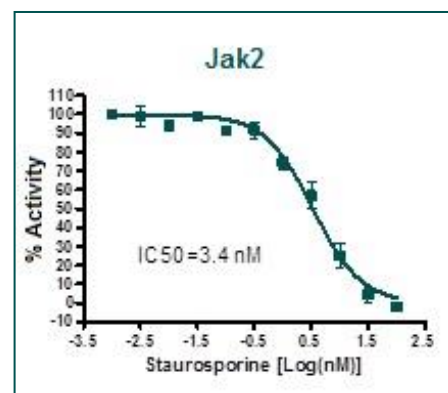
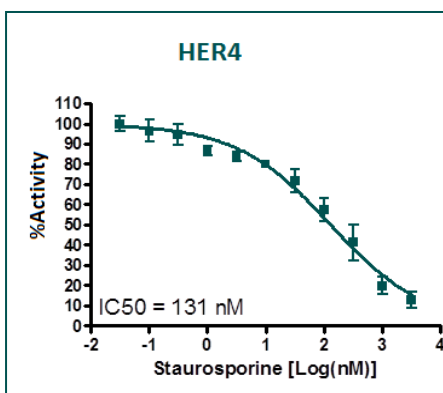
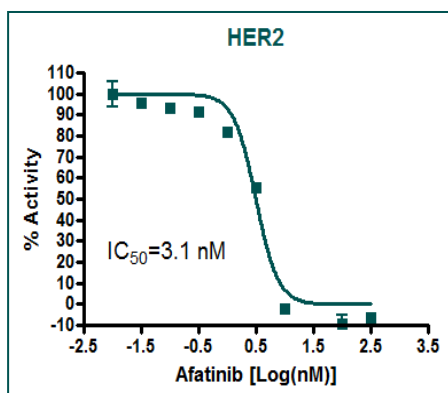
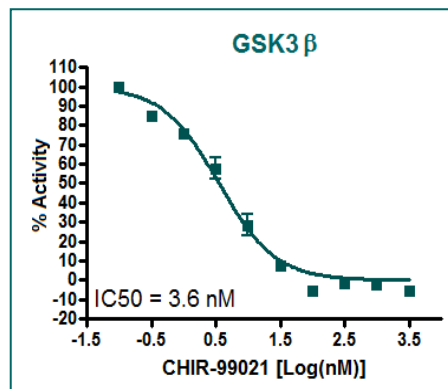
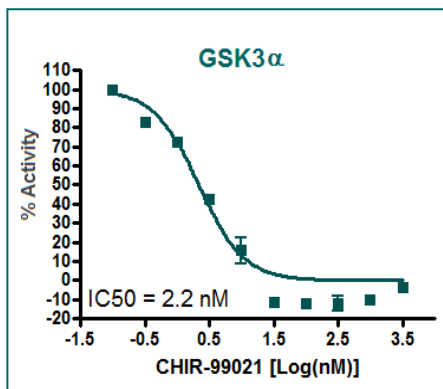
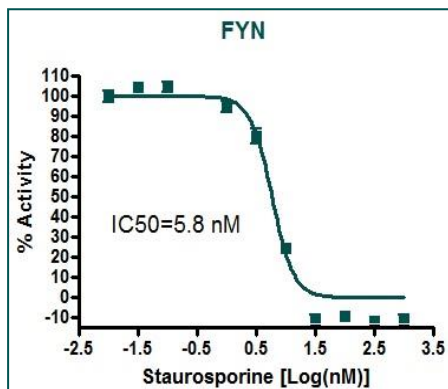
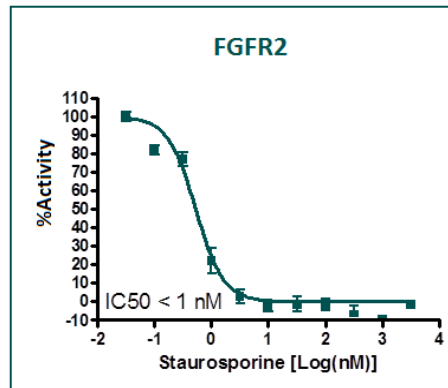
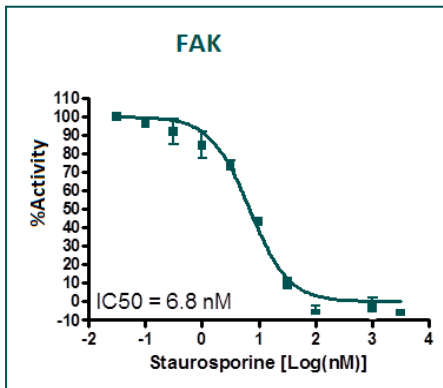
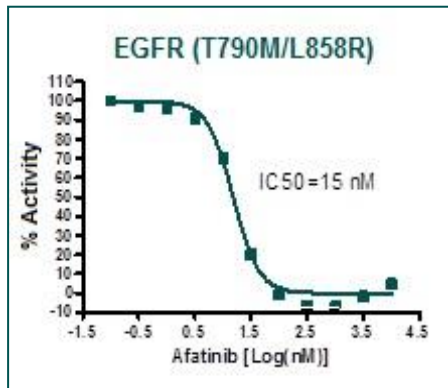


OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

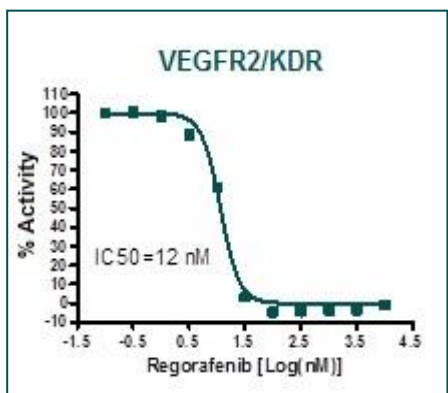
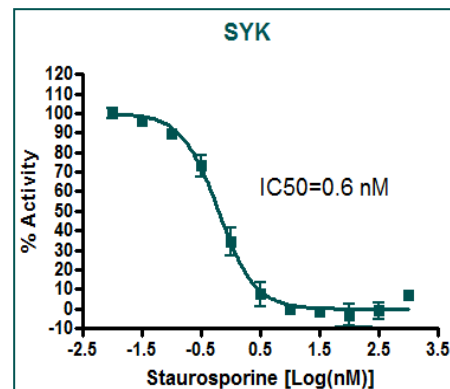
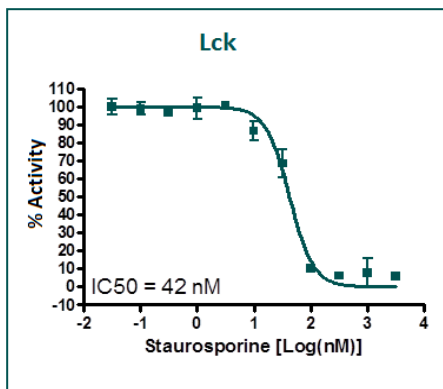
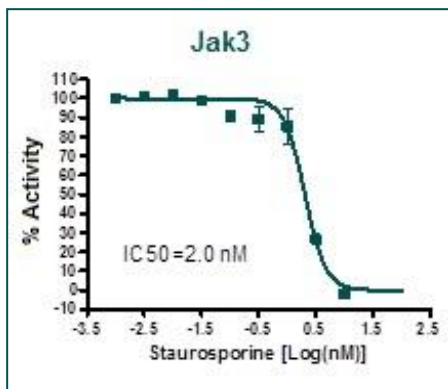


OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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



Our panel of kinases includes:

ALK	FAK
BRAF	FGFR2
BRAF (V600E)	FYN
BTK	GSK3 α
c-Src	GSK3 β
CDK4/CyclinD3	Her2
EGFR (WT)	Her4
EGFR (L858R)	Lck
EGFR (T790M)	Jak2
EGFR (L858R, T790M)	Jak3
EGFR (murine)	SYK
	VEGFR2/KDR

Application Reference(s): Figueroa, M., Oberlies, N.H., *et al.*, *The Journal of Antibiotics*. **65**, 559-564 (2012).

Also available: **Metabolic Kinase Profiling!** Pyruvate Kinases (PKM1, PKM2, PKR, PKL), Fucokinase, Galactokinase, and more. Contact us for details.

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