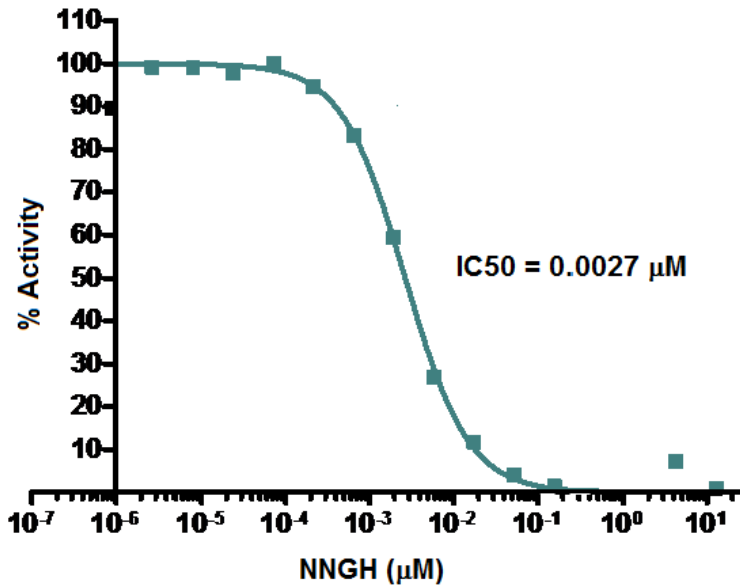
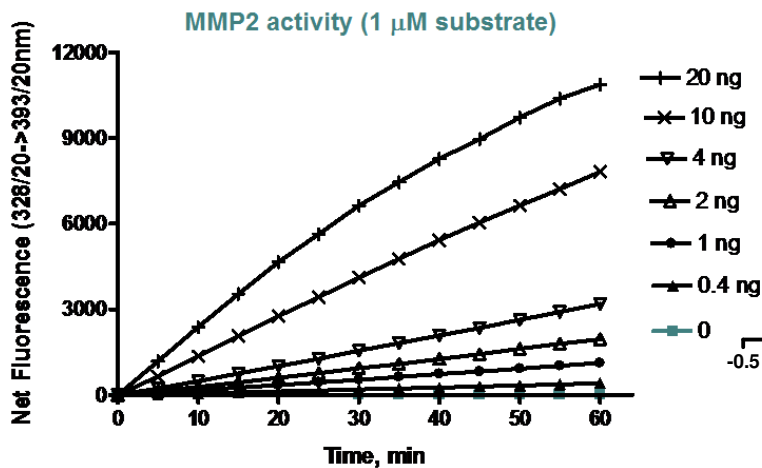


MMP2 Screening Services

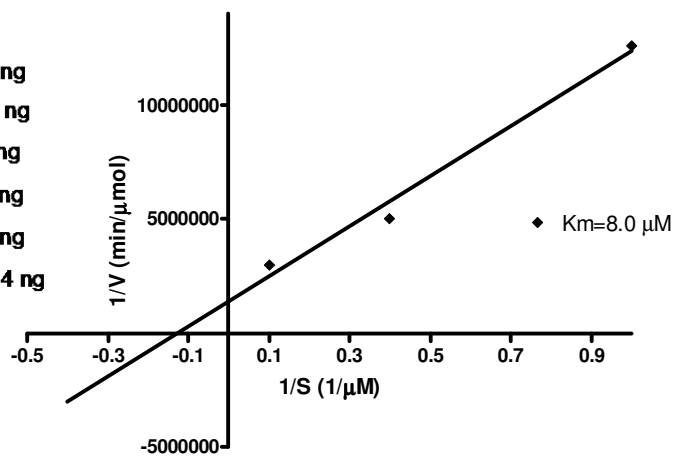


Please contact us for more information on our screening and profiling services.

Kinetics studies



MMP2 Lineweaver-Burk plot



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