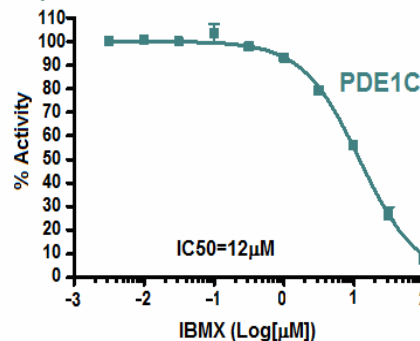
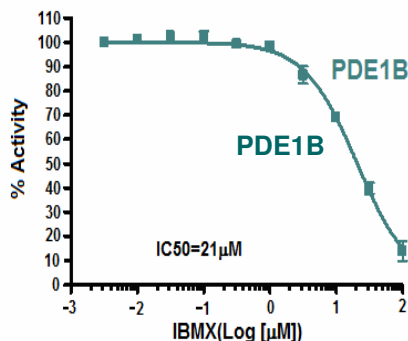
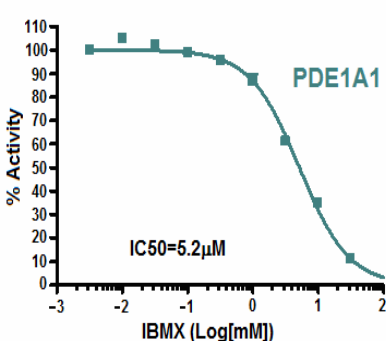


PDE Profiling

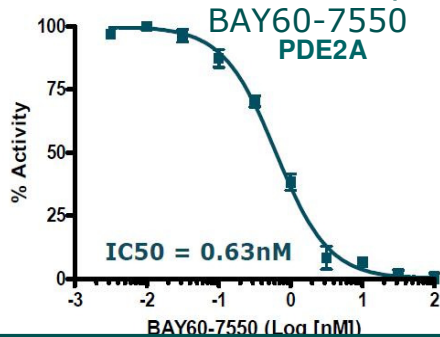
Over 22 different PDE isozymes available!

Contact us for more information on our screening and profiling services.

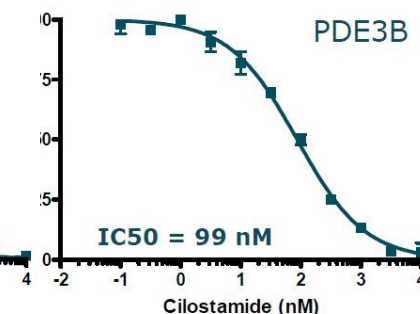
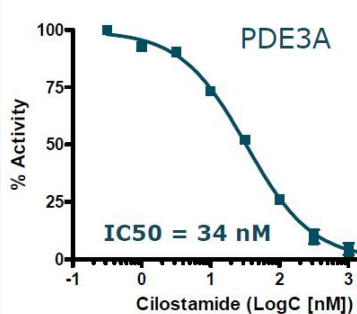
PDE1A, PDE1B, and PDE1C inhibition by IBMX



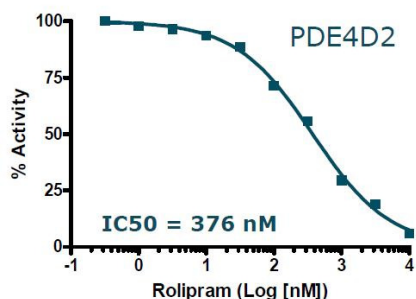
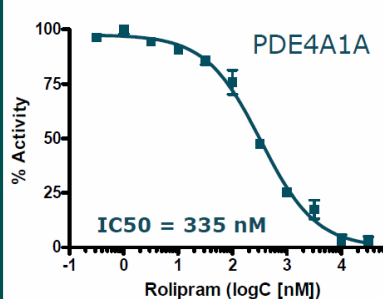
PDE2A inhibition by BAY60-7550



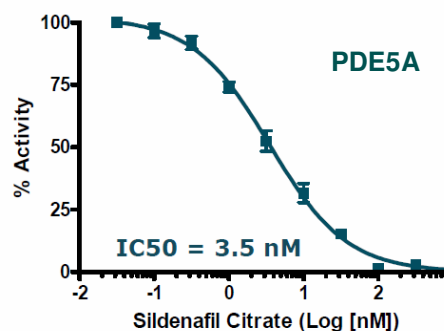
PDE3A and PDE3B inhibition by Cilostamide



PDE4A and PDE4D inhibition by Rolipram



PDE5A inhibition by Sildenafil



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