



6405 Mira Mesa Blvd Ste 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## **Data Sheet**

### ***Terbium (Tb) Donor*** **Catalog #60394**

**DESCRIPTION:** The Terbium Donor, also known as Tb Donor, is used for phosphodiesterase (PDE) TR-FRET (Time Resolved Fluorescence Resonance Energy Transfer) technology based assays including PDE1B, PDE1C, PDE3A, PDE4D, and PDE10A. In these assays, terbium conjugates with phosphate.

**PURITY:** >95% by TLC.

**APPLICATIONS:** Suitable for studying phosphodiesterase activity and for screening small molecule inhibitors for drug discovery and HTS applications.

**SUPPLIED AS:** An aqueous solution

**QUANTITY:** 50  $\mu$ l

**STORAGE:** -20°C

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)