



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@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: 30 μ l @ 2 μ M

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 **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com