

R&D, Administration and Europe Office  
Cisbio Bioassays  
Phone: +33 (0)4 66 79 67 05  
Fax: +33 (0)4 66 79 19 20  
E-mail: bioassays@cisbio.com

Japan Office  
Sceti Medical Labo K.K.  
Phone: +81 (0)3 5510 2932  
Fax: +81 (0)3 5510 0130  
E-mail: reagent@scetimedilabo.co.jp

USA Office  
Cisbio US, Inc.  
Phone : 888 963 4567  
Fax : 781 687 1500  
E-mail : htrfi nfo@cisbio.us

[www.htrf.com](http://www.htrf.com)

## Product Information

For in vitro research use only  
Version June 2010 - A



### Tag-lite® CLIP Lumi4® -Tb

<b>Product number:</b>	<b>SCLPTBE</b>	<b>20 nmoles</b>
	<b>SCLPTBF</b>	<b>5 x 20 nmoles</b>
	<b>SCLPTBZ</b>	<b>500 nmoles</b>

Storage temperature: -20°C

**Formulation** Dessicated form

## Description

This substrate is a CLIP Substrate labeled with a Lumi4-Tb.

## Reconstitution

Vial of 20 nmoles : add 200 µl DMSO to get a 100 µM solution.  
Vial of 500 nmoles : add 5 ml DMSO to get a 100 µM solution.

## Storage and stability

Upon receiving the ligand, store at -20°C or colder .  
After reconstitution with DMSO, product must be used immediately or dispensed into aliquots and frozen at -80°C. Do not repeat freezing and thawing.

## Protocol

Information available on: [www.htrf.com](http://www.htrf.com)

## MSDS

Information available on: [www.htrf.com](http://www.htrf.com)