

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

Product Information

For in vitro research use only
Version June 2010 - A

Tag-lite® SNAP Lumi4® -Tb

Product number:	SSNPTBC	2 nmoles
	SSNPTBD	5 nmoles
	SSNPTBG	5 x 5 nmoles
	SSNPTBX	100 nmoles

Storage temperature: -20°C

Formulation Dessicated form

Description

This substrate is a SNAP Substrate labeled with a Lumi4-Tb

Reconstitution

Vial of 2 nmoles : add 20 µl DMSO to get a 100 µM solution
Vial of 5 nmoles : add 50 µl DMSO to get a 100 µM solution
Vial of 100 nmoles : add 1 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

MSDS

Information available on: www.htrf.com