

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 Red

Product number:	SSNPREDE	20 nmoles
	SSNPREDF	5 x 20 nmoles
	SSNPREDZ	500 nmoles

Storage temperature: -20°C

Formulation Dessicated form

Description

This substrate is a SNAP Substrate labeled with a red emitting HTRF fluorescent probe.

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

MSDS

Information available on: www.htrf.com