



Tag-lite® SNAP-green

Product Information: #2 September 2023

- B

Part #: SSNPGRNE 5 nmoles
SSNPGRNZ 100 nmoles

Store at: -20°C

Formulation: Dessicated form

For research use only. Not for use in diagnostic procedures.

DESCRIPTION

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

RECONSTITUTION

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 substrate, 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.

MSDS

Information available on: www.revvity.com



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

www.revvity.com

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.