



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data Sheet

PL-01 Assay Buffer

Catalog #: 33298

DESCRIPTION: The *PL-01 Assay Buffer* is optimized for use with the following BPS Bioscience assay kits:

- | | | |
|--|--------|---------------|
| • PCSK9[Biotinylated]-LDLR Binding Assay Kit | #72002 | 96 reactions |
| • PCSK9-LDLR TR-FRET Assay Kit | #72010 | 384 reactions |
| • PCSK9(D374T)-LDLR TR-FRET Assay Kit | #72011 | 384 reactions |

APPLICATIONS: Studying enzyme kinetics and PCSK9 applications in drug discovery.

QUANTITY: 50 ml

STORAGE: -20°C. Avoid freeze/thaw cycles.

STABILITY: Stable for at least one year from date of receipt when stored as directed.

INSTRUCTIONS FOR USE: For a detailed protocol on the use of this product, please see the protocol for PCSK9-LDLR TR-FRET Assay Kit (BPS Bioscience #72010).

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
PCSK9, His-tag	71304	20 µg
PCSK9, Biotin	71211	20 µg
Anti-PCSK9 Antibody	71207	50 µg
LDLR, FLAG-tag	71205	50 µg
PCSK9-LDLR TR-FRET Kit	72010	384 rxns
PCSK9(D374T)-LDLR TR-FRET Kit	72011	384 rxns

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com