

Product Information



TR-FRET Assay Buffer 2 (10X)

Item No. 600508

IMPORTANT

Water used to prepare all reagents and buffers must be deionized and free of trace organic contaminants ('UltraPure'). Use activated carbon filter cartridges or other organic scavengers. Glass distilled water (even if double distilled), HPLC-grade water, and sterile water (for injections) are not adequate for TR-FRET. UltraPure water may also be purchased (Item No. 400000).

Laboratory Procedures

2 ml TR-FRET Assay Buffer 2 (10X):

This vial contains 2 ml TR-FRET Assay Buffer 2 (10X). This assay buffer, when stored as supplied, is stable for one year at -20°C. Prior to use, add 18 ml of UltraPure water to the vial. For best results, filter completed 1X assay buffer with a 0.22 µm filter before use. *NOTE: It is normal for the concentrated buffer to contain crystalline salts after thawing. These will completely dissolve upon dilution with water.* Store the buffer at 4°C; it will be stable for approximately one month.

OR

10 ml TR-FRET Assay Buffer 2 (10X):

This vial contains 10 ml TR-FRET Assay Buffer 2 (10X). This assay buffer, when stored as supplied, is stable for one year at -20°C. Prior to use, dilute to a total volume of 100 ml with UltraPure water. For best results, filter completed 1X assay buffer with a 0.22 µm filter before use. *NOTE: It is normal for the concentrated buffer to contain crystalline salts after thawing. These will completely dissolve upon dilution with water.* Store the buffer at 4°C; it will be stable for approximately one month.

OR

50 ml TR-FRET Assay Buffer 2 (10X):

This vial contains 50 ml TR-FRET Assay Buffer 2 (10X). This assay buffer, when stored as supplied, is stable for one year at -20°C. Prior to use, dilute to a total volume of 500 ml with UltraPure water. For best results, filter completed 1X assay buffer with a 0.22 µm filter before use. *NOTE: It is normal for the concentrated buffer to contain crystalline salts after thawing. These will completely dissolve upon dilution with water.* Store the buffer at 4°C; it will be stable for approximately one month.

CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with the L3MBTL1 MBT Domains TR-FRET Assay Kit (Item No. 601030) from Cayman. This product may not perform as described if used with other assay reagents or protocols.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/600508

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/27/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com