TACE substrate II (fluorogenic)

Sensitive fluorogenic substrate for TACE (TNF α converting enzyme, ADAM17). Also cleaved by ADAM10 and likely other ADAMs and MMPs. The k_{cat}/K_{m} of TACE for this substrate is 7-fold higher than for the same peptide with the Mca-Dpa fluorophore-quencher pair. FAM fluorescence is quenched by the TAMRA group until cleavage separates them. Ex.: 485nm, Em.: 535nm, although the following Ex/Em can also be used: 490,494/515,520nm. This substrate is useful for inhibitor screening, kinetic analysis, and cellular activity assay.

Citations: 1

View Online »

Ordering Information

Order Online »

BML-P228-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light and moisture.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ADAM17 substrate II (fluorogenic)

Appearance Solid.

Formulation Lyophilized.

MW 2157.3

Purity ≥95% (HPLC)

Sequence 5-FAM-Ser-Pro-Leu-Ala-Gln-Ala-Val-Arg-Ser-Ser-Ser-Arg-

Lys(5-TAMRA)-NH₂

Solubility Soluble in DMSO.

Last modified: May 29, 2024

