

PRODUCT INFORMATION



Dabcyl-K-SARS-CoV-2 Replicase Polyprotein 1ab(3235-3246)-E-EDANS (trifluoroacetate salt)

Item No. 37409

Formal Name: N²-[4-[[4-(dimethylamino)phenyl]azo]benzoyl]-L-lysyl-L-threonyl-L-seryl-L-alanyl-L-valyl-L-leucyl-L-glutamyl-L-serylglycyl-L-phenylalanyl-L-arginyl-L-lysyl-L-methionyl-N-[2-[[5-sulfo-1-naphthalenyl]amino]ethyl]-L-α-glutamine, trifluoroacetate salt

Synonyms: Dabcyl-KTSAVLQSGFRKME-EDANS, Dabcyl-Lys-Thr-Ser-Ala-Val-Leu-Gln-Ser-Gly-Phe-Arg-Lys-Met-Glu-EDANS

MF: C₉₅H₁₄₁N₂₅O₂₄S₂ • XCF₃COOH

FW: 2,081.4

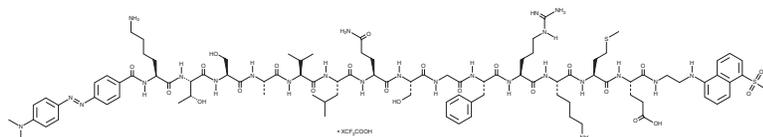
Purity: ≥90%

Ex./Em. Max: 340/485 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Dabcyl-K-SARS-CoV-2 replicase polyprotein 1ab (3235-3246)-E-EDANS (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Dabcyl-K-SARS-CoV-2 replicase polyprotein 1ab (3235-3246)-E-EDANS (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Dabcyl-K-SARS-CoV-2 replicase polyprotein 1ab (3235-3246)-E-EDANS (trifluoroacetate salt) is soluble in DMSO.

Description

Dabcyl-K-SARS-CoV-2 replicase polyprotein 1ab (pp1ab) (3235-3246)-E-EDANS is a fluorogenic substrate for severe acute respiratory syndrome coronavirus (SARS-CoV) and SARS-CoV-2 main protease (M^{pro}), also known as 3C-like protease (3CL^{pro}).^{1,2} Upon enzymatic cleavage by SARS-CoV or SARS-CoV-2 M^{pro} at the Gln-Ser site, EDANS is separated from the Dabcyl quencher and its fluorescence can be used to quantify SARS-CoV or SARS-CoV-2 M^{pro} activity.¹⁻³ EDANS displays excitation/emission maxima of 340/485 nm, respectively.

References

1. Hsu, J.T.-A., Kuo, C.-J., Hsieh, H.-P., *et al.* Evaluation of metal-conjugated compounds as inhibitors of 3CL protease of SARS-CoV. *FEBS Lett.* **574(1-3)**, 116-120 (2004).
2. Xiao, T., Wei, Y., Cui, M., *et al.* Effect of dihydromyricetin on SARS-CoV-2 viral replication and pulmonary inflammation and fibrosis. *Phytomedicine* **91**, 153704 (2021).
3. Kim, Y., Lovell, S., Tiew, K.-C., *et al.* Broad-spectrum antivirals against 3C or 3C-like proteases of picornaviruses, noroviruses, and coronaviruses. *J. Virol.* **86(21)**, 11754-11762 (2012).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/09/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM