

PRODUCT INFORMATION



MDM2 (32-46) (trifluoroacetate salt)

Item No. 38836

Formal Name: ((S)-2-formylpyrrolidin-1-yl)-L-leucyl-L-leucyl-L-leucyl-L-lysyl-L-leucyl-L-leucyl-L-leucyl-L-lysyl-L-seryl-L-valylglycyl-L-alanyl-L-glutaminy-L-lysyl-L-aspartic acid, trifluoroacetate salt

Peptide Sequence: PLLLLKLLKSVGAQKD-OH

MF: $C_{75}H_{135}N_{19}O_{20} \bullet XCF_3COOH$

FW: 1,623.0

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

H—Pro—Leu—Leu—Leu—Lys—Leu—Leu—Lys—Ser—Val—
Gly—Ala—Gln—Lys—Asp—OH
 $\bullet XCF_3COOH$

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

Laboratory Procedures

MDM2 (32-46) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the MDM2 (32-46) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. MDM2 (32-46) (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of MDM2 (32-46) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of MDM2 (32-46) (trifluoroacetate salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

MDM2 (32-46) is a peptide fragment of MDM2.

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, 04/24/2024

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