

PRODUCT INFORMATION



HOXB7 (8-25) (human)

Item No. 41334

Formal Name: (S)-N-(((S)-1-(((S)-6-amino-1-(((S)-1-(((S)-1-(((2R,3S,6S)-6,8-diamino-2-hydroxy-4,5,8-trioxooctan-3-yl)amino)-4-methyl-1-oxopentan-2-yl)amino)-1-oxo-3-phenylpropan-2-yl)amino)-3-hydroxy-1-oxopropan-2-yl)amino)-1-oxohexan-2-yl)amino)-3-(4-hydroxyphenyl)-1-oxopropan-2-yl)-1-(hydroxy-L-alanylglycyl-L-threonyl-L-alanyl-L-phenylalanyl-L-valyl-L-seryl-L-seryl-L-seryl-L-alanyl)pyrrolidine-2-carboxamide

H—Asn—Thr—Leu—Phe—Ser—Lys—Tyr—Pro—Ala—Ser—
Ser—Ser—Val—Phe—Ala—Thr—Gly—Ala—OH

Synonym: Homeobox Protein B7 (8-25)

Peptide Sequence: NTLFSKYPASSVFATGA-OH

MF: $C_{84}H_{126}N_{20}O_{27}$

FW: 1,848

Purity: ≥95%

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

HOXB7 (8-25) (human) is supplied as a solid. A stock solution may be made by dissolving the HOXB7 (8-25) (human) in the solvent of choice, which should be purged with an inert gas. HOXB7 (8-25) (human) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

HOXB7 (8-25) (human) is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

HOXB7 (8-25) is a peptide fragment of the transcription factor homeobox protein B7 (HOXB7).

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, 07/24/2024

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