

PRODUCT INFORMATION

Histone H2B (21-35) (human, mouse, rat, bovine) (trifluoroacetate salt)
Item No. 27472

Formal Name: L-alanyl-L-glutamyl-L-lysyl-L-lysyl-L-aspartylglycyl-L-lysyl-L-lysyl-L-arginyl-L-lysyl-L-arginyl-L-seryl-L-arginyl-L-lysyl-L-glutamic acid, trifluoroacetate salt

Synonyms: AQKKDGKKRKRSRKE, AspH2B₂₁₋₃₅, H—Ala—Gln—Lys—Lys—Asp—Gly—Lys—Lys—Arg—Lys—Arg—Ser—Arg—Lys—Glu—OH
H2B₂₁₋₃₅ • XCF₃COOH

MF: C₇₆H₁₄₃N₃₁O₂₂ • XCF₃COOH

FW: 1,843.2

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

Histone H2B (21-35) (human, mouse, rat, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H2B (21-35) (human, mouse, rat, bovine) (trifluoroacetate salt) in water. The solubility of histone H2B (21-35) (human, mouse, rat, bovine) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histone H2B (21-35) is a peptide fragment of histone H2B that corresponds to amino acid residues 22-36 of the human histone H2B sequence. Histone H2B (21-35) contains an aspartic acid at position 25 that spontaneously converts to isoaspartic acid at physiological pH *in vitro* and *in vivo* and induces production of autoantibodies against histone H2B in murine models of systemic lupus erythematosus (SLE).¹

Reference

1. Doyle, H.A., Aswad, D.W., and Mamula, M.J. Autoimmunity to isomerized histone H2B in systemic lupus erythematosus. *Autoimmunity* **46**(1), 6-13 (2013).

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, 11/30/2022

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