

PRODUCT INFORMATION

Histone H2AX (134-143) (human) (trifluoroacetate salt)

Item No. 27441

Formal Name:	L-lysyl-L-lysyl-L-alanyl-L-threonyl-L-glutaminyl-L-alanyl-L-seryl-L-glutaminyl-L- α -glutamyl-L-tyrosine, trifluoroacetate salt
Synonyms:	H-Lys-Lys-Ala-Thr-Gln-Ala-Ser-Gln-Glu-Tyr-OH, KKATQASQEY
MF:	$C_{49}H_{80}N_{14}O_{18} \cdot XCF_3COOH$
FW:	1,153.3
Purity:	$\geq 95\%$
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 H2AX (134-143) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H2AX (134-143) (human) (trifluoroacetate salt) in water. The solubility of histone H2AX (134-143) (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histone H2AX (134-143) is a C-terminal peptide fragment of histone H2AX that corresponds to amino acid residues 134-143 of the human histone H2AX sequence. Histone H2AX (134-143) is an evolutionarily conserved motif that differentiates histone H2AX from other H2A histones.¹ It contains a serine residue that is rapidly phosphorylated upon introduction of DNA double-strand breaks and initiation of DNA fragmentation during apoptosis.

Reference

1. Rogakou, E.P., Nieves-Neira, W., Boon, C., *et al.* Initiation of DNA fragmentation during apoptosis induces phosphorylation of H2AX histone at serine 139. *J. Biol. Chem.* **275**(13), 9390-9395 (2000).

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