

PRODUCT INFORMATION



TAF 10 Peptide

Item No. 10228

Synonyms: TAF10 RNA Polymerase II, TATA Box Binding Protein (TBP)-Associated Factor, TBP-Associated Factor 10, TAF_{II}30

Peptide Sequence: Human TAF10 amino acids 186-195; (Ac)SKSKDRKYTL

FW: 1,267

Stability: ≥3 years at -20°C

Supplied as: A 1 mg peptide lyophilized from bicarbonate buffer

Solubility: ≥20 mg/ml with water or aqueous buffers

Description

TAF10 is one of many protein factors or coactivators associated with RNA polymerase II activity.¹ One vial of this peptide may be used as a methyltransferase acceptor peptide for more than 200 reactions at 15 μM.^{2,3}

References

1. Metzger, D., Scheer, E., Soldatov, A., *et al.* Mammalian TAF_{II}30 is required for cell cycle progression and specific cellular differentiation programmes. *EMBO J.* **18(17)**, 4823-4834 (1999).
2. Couture, J.-F., Collazo, E., Hauk, G., *et al.* Structural basis for the methylation site specificity of SET7/9. *Nature Structural and Molecular Biology* **13(2)**, 140-146 (2006).
3. Collazo, E., Couture, J.-F., Bulfer, S., *et al.* A coupled fluorescent assay for histone methyltransferases. *Anal. Biochem.* **342**, 86-92 (2005).

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, 11/06/2023

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