

Product Information



TAF10 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:	≥1 year at -20°C
Supplied as:	1 mg peptide lyophilized from bicarbonate buffer
Solubility:	At least 20 mg/ml with water or aqueous buffers

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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10228

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 04/09/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com