

Product Information



EAF2 Blocking Peptide

Catalog No. 10005732

EAF2 is a testosterone-regulated apoptosis inducer with tumor suppressive activity.¹ *In vivo*, overexpression of EAF2 induces massive apoptosis and inhibits prostate tumor growth.¹ It was first identified as ELL-associated factor 2, which shares high homology with EAF1 that also interacts with ELL, and may play a role in leukemogenesis.² EAF2 mRNA is widely expressed in various tissues, with the most abundant expression in the prostate, kidney, bone marrow, and lymph nodes.¹ Cayman's EAF2 polyclonal antibody detects a 43 kDa endogenous protein by western blot analysis in human, murine, and rat samples, which is consistent with published data for EAF2.²

Laboratory Procedures

This vial contains 200 µg peptide in 200 µl TBS, pH 7.4, containing 0.1% BSA and 0.02% sodium azide. The peptide is derived from the human EAF2 sequence amino acids 5-16 (AGFSLDRRERV). Store aliquots at -20°C; they will be good for up to two years. This blocking peptide can be used in conjunction with Cayman's EAF2 Polyclonal Antibody (Item No. 10005190) to block protein-antibody complex formation during immunochemical analysis of EAF2.

Reconstitute the lyophilized peptide with 200 µl of PBS or distilled water. Store this peptide solution at -20°C. It will be stable for at least two years. To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the EAF2 Polyclonal Antibody (Item No. 10005190) and blocking peptide together in a 1:1 (v/v) ratio in a microfuge tube. For example, mix 20 µl of antibody and 20 µl of peptide.*
2. Incubate for 1 hour at room temperature with occasional mixing prior to further dilution and application of the mixture to the immunoblot.
3. Dilute the mixture to the final working antibody concentration and apply to the slide or membrane as usual.

*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur.

References

1. Xiao, W., Zhang, Q., Jiang, F., *et al.* Suppression of prostate tumor growth by U19, a novel testosterone-regulated apoptosis inducer. *Cancer Res.* **63**, 4698-4704 (2003).
2. Simone, F., Luo, R.T., Polak, P.E., *et al.* ELL-associated factor 2 (EAF2), a functional homolog of EAF1 with alternative ELL binding properties. *Blood* **101**(6), 2355-2362 (2003).

Related Product

EAF2 Polyclonal Antibody - Cat. No. 10005190

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

MATERIAL 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 Material 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, 10/05/2010

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