

Product Information



EAF2 Polyclonal Antibody

Item No. 10005190

- Synonyms:** ELL-Associated Factor 2, U19
- Contents:** This vial contains 150 µg peptide affinity-purified IgG in 300 µl TBS, pH 7.4, containing 50% glycerol, 0.1% BSA, and 0.02% sodium azide
- Host:** Rabbit
- Antigen:** Human EAF2 amino acids 5-16. The antigen alignment with other sequences is as follows:
Human AGFSHLDRRERV
Mouse AG1a y LDRRERV
Rat AG1a y LDRRER i
- Cross Reactivity:** (+) Human, mouse, and rat EAF2; other species not tested
- Stability:** ≥1 year at -20°C
- Applications:** Recommended starting dilutions: western blot; 1.5 µg/ml and immunohistochemistry (formalin-fixed paraffin-embedded sections); 2.5 µg/ml. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

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, mouse, and rat samples, which is consistent with published data for EAF2.²

References

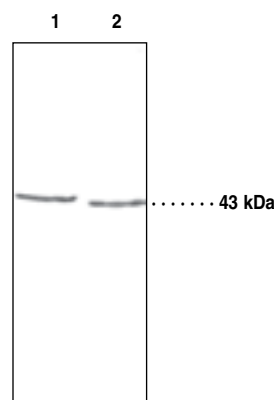
- Xiao, W., Zhang, Q., Jiang, F., *et al.* Suppression of prostate tumor growth by U19, a novel testosterone-regulated apoptosis inducer. *Cancer Res.* **63**, 4698-4704 (2003).
- Simone, F., Luo, R.T., Polak, P.E., *et al.* ELL-associated factor 2 (EAF2), a functional homolog of EAF1 with alternative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

Laboratory Procedures

Immunoperoxidase immunohistochemical procedure

A. Paraffin sections

- Deparaffinize sections 3 times with xylene or a xylene substitute, 5 minutes each.
- Rehydrate sections 2 times with 100% ethanol, 5 minutes each, followed by 95%, 90%, 80%, 70%, 50% ethanol, 5 minutes each.
- Rinse sections in distilled water for 5 minutes.
- Block endogenous peroxidase activity with 0.3% H₂O₂ in water (use methanol instead of water in case of strong endogenous peroxidase activity) for 15 minutes.
- Wash sections 3 times in TBS, pH 7.4, 5 minutes each.
- Incubate sections with 5% normal serum (same species as secondary antibody host) for 30 minutes.
- Rinse sections 3 times with TBS, pH 7.4, 5 minutes each.
- Incubate sections with 2.5 µg/ml polyclonal antibody (recommended starting dilution; optimal dilution to be determined by end user) for 1 hour at room temperature (in a humid chamber if desired).
- Wash sections 3 times in TBS, pH 7.4, 5 minutes each.
- Incubate sections for 30 minutes with biotinylated secondary antibody using a dilution as recommended by provider.
- Wash sections 3 times in TBS, pH 7.4, 5 minutes each.
- Incubate sections for 30 minutes with ABC reagent using a dilution as recommended by provider.



Lane 1: Rat liver 100 k x g pellet re-suspension (50 µg)
Lane 2: Murine liver 10 k x g pellet re-suspension (100 µg)

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, 12/20/2011

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



13. Wash sections 3 times in TBS, pH 7.4, 5 minutes each.
14. Incubate sections in peroxidase substrate solution. Check staining under a microscope frequently. When desired staining intensity is achieved, rinse sections with distilled water thoroughly.
15. Counter stain sections if desired. Rinse sections thoroughly after counter stain.
16. Dehydrate sections through 50%, 70%, 80%, 90%, 95%, 100% (2 times) ethanol for 5 minutes each.
17. Clear sections 3 times with xylene or xylene substitute, 5 minutes.
18. Mount sections with coverslips.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10005190

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