

# Product Information

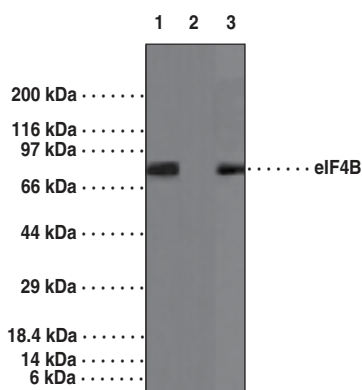


## eIF4B Peptide Affinity-Purified Polyclonal Antibody (aa 570-630)

Item No. 13529

- Contents:** This vial contains 100 µg of peptide affinity-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
- Synonym:** Eukaryotic Translation Initiation Factor-4B
- Antigen:** Synthetic peptide from human eIF4B within the region of amino acids 570-630
- Host:** Rabbit
- Cross Reactivity:** (+) Chimpanzee, human, mouse, orangutan, and Rhesus monkey eIF4B
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Applications:** The recommended starting dilution for western blot is 0.005-1.0 µg/ml.

Eukaryotic initiation factor-4B peptide (eIF4B) is a multidomain eIF with a range of activities. It contains one RNA recognition motif (RRM) domain, a consensus RNA-binding site (RNP-CS) near the N terminus, and a central domain rich in aspartic acid, arginine, tyrosine, and glycine (DRYG). The DRYG domain of eIF4B helps in self-association and interaction with the p170 subunit of eIF3. eIF4B binds to both mRNA and 18S ribosomal RNA and thus acts as a bridge between the mRNA and the 40S ribosomal subunit during translation initiation. The ATPase and RNA helicase activity of the DEAD box protein eIF4A is stimulated by eIF4B and thus functions in close association with eIF4E, a heterotrimeric protein that binds to mRNA cap structure and unwinds inhibitory secondary structure in the 5' untranslated region of eucaryotic mRNAs in an ATP-dependent manner. Ubiquitous expression is seen in most tissues.



Lane 1: K562 cell lysate (Absence of immunizing peptide)  
Lane 2: K562 cell lysate (Presence of immunizing peptide)  
Lane 3: RAW cell lysate

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/13529](http://www.caymanchem.com/catalog/13529)

### 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](http://www.caymanchem.com)

**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, 12/15/2011