

Product Information

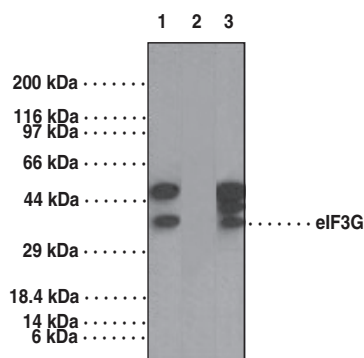


eIF3G Polyclonal Antibody (aa 200-250)

Catalog No. 13537

- 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-3G
- Antigen:** Synthetic peptide from human eIF3G within the region of amino acids 200-250
- Host:** Rabbit
- Cross Reactivity:** (+) Chimpanzee, bovine, canine, equine, human, murine, rat, and Rhesus monkey eIF3G
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Applications:** The recommended starting dilution for western blot is 0.025-0.1 µg/ml and 5 µg/ml for immunohistochemistry (paraffin-embedded sections).

Eukaryotic translation initiation factor-3G (eIF3G) is an RNA-associated protein involved in protein synthesis. It is a member of the eIF3 subunit G family and the largest of the eIF family with eIF3 being a multiprotein complex composed of at least 13 different subunits. eIF3G has a RNA recognition motif (RRM) domain and is known to interact with eIF3K. The function of eIF3G is to bind to the 40S ribosomal subunit and promote the binding of methionyl-tRNAi and mRNA. This subunit of eIF3 also binds to the 18S rRNA. It is ubiquitously expressed in most tissues.



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

Related Products

eIF4B Polyclonal Antibody (aa 570-630) - Cat. No. 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

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, 07/14/2010