

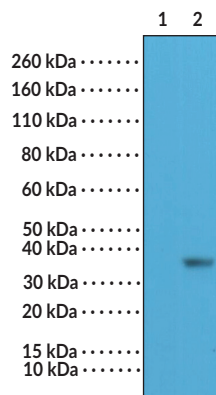
# PRODUCT INFORMATION

**eIF2 $\alpha$  (Phospho-Ser<sup>51</sup>) Rabbit Monoclonal Antibody (Clone RM298)**  
 Item No. 32244

## Overview and Properties

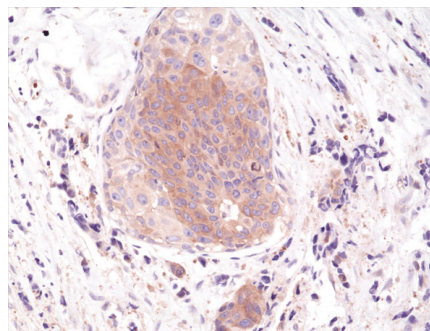
**Contents:** This vial contains 100  $\mu$ l of protein A-affinity purified monoclonal antibody.  
**Synonyms:** Eukaryotic Translation Initiation Factor 2 Subunit 1  
**Immunogen:** Phosphopeptide corresponding to human eIF2 $\alpha$  (phospho-Ser<sup>51</sup>)  
**Cross Reactivity:** (+) eIF2 $\alpha$ ; (-) eIF2 $\alpha$  without phosphorylation at serine 51  
**Species Reactivity:** (+) Human  
**Form:** Liquid  
**Storage:** -20°C (as supplied)  
**Stability:**  $\geq$ 1 year  
**Storage Buffer:** PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide  
**Clone:** RM298  
**Host:** Rabbit  
**Isotype:** IgG  
**Applications:** Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

## Images

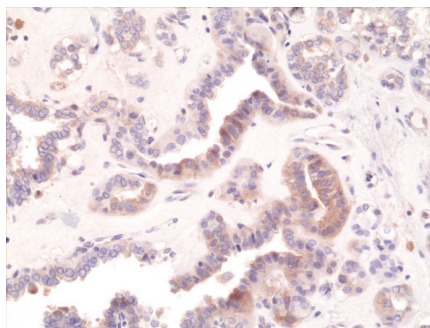


**Lane 1:** HeLa cell lysates untreated  
**Lane 2:** HeLa cell lysates treated

WB of HeLa cell lysates untreated or treated with calyculin A (Item No. 19246) using eIF2 $\alpha$  (Phospho-Ser<sup>51</sup>) Rabbit Monoclonal Antibody (Clone RM298) at a dilution of 1:200.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast cancer tissue using eIF2 $\alpha$  (Phospho-Ser<sup>51</sup>) Rabbit Monoclonal Antibody (Clone RM298) at a 1:200 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human thyroid cancer tissue using eIF2 $\alpha$  (Phospho-Ser<sup>51</sup>) Rabbit Monoclonal Antibody (Clone RM298) at a 1:200 dilution.

**WARNING**  
 THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
 This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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**  
 Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/30/2024

**CAYMAN CHEMICAL**  
 1180 EAST ELLSWORTH RD  
 ANN ARBOR, MI 48108 · USA  
 PHONE: [800] 364-9897  
 [734] 971-3335  
 FAX: [734] 971-3640  
 CUSTSERV@CAYMANCHEM.COM  
 WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

eIF2 $\alpha$  is the  $\alpha$  subunit of eukaryotic inhibition factor 2, which also includes eIF2 $\beta$  and eIF2 $\gamma$  subunits and is involved in the initiation of translation.<sup>1</sup> eIF2 $\alpha$  is composed of a globular N-terminal domain, which contains an S1 domain and an  $\alpha$ -helical domain, and a globular C-terminal domain.<sup>2</sup> eIF2 $\alpha$  can be phosphorylated at serine 51 by the eIF2 $\alpha$  kinases HRI/EIF2AK1, PKR/EIF2AK2, PERK/EIF2AK3, and GCN2/EIF2AK4 following various cell and ER stressors, viral invasion, or nutrient deprivation, among other factors.<sup>3-5</sup> Phosphorylation of eIF2 $\alpha$  at serine 51 promotes the interaction between eIF2 $\alpha$  and the guanine nucleotide exchange factor eIF2B, which inhibits eIF2B activity and prevents the initiation of translation, reducing global protein synthesis in the cell.<sup>1</sup> The transcription of specific factors involved in the integrated stress response (ISR), such as ATF4, is upregulated by eIF2 $\alpha$  phosphorylation.<sup>3</sup> Dephosphorylation of eIF2 $\alpha$  by the protein phosphatase 1 (PP1) complex, which includes GADD34 and CREP, terminates the ISR and returns protein synthesis back to basal levels. Cayman's eIF2 $\alpha$  (Phospho-Ser<sup>51</sup>) Rabbit Monoclonal Antibody (Clone RM298) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

## References

---

1. Sudhakar, A., Ramachandran, A., Ghosh, S., *et al.* Phosphorylation of serine 51 in initiation factor 2 $\alpha$  (eIF2 $\alpha$ ) promotes complex formation between eIF2 $\alpha$ (P) and eIF2B and causes inhibition in the guanine nucleotide exchange activity of eIF2B. *Biochemistry* **39**(42), 12929-12938 (2000).
2. Ito, T., Marintchev, A., and Wagner, G. Solution structure of human initiation factor eIF2 $\alpha$  reveals homology to the elongation factor eEF1B. *Structure* **12**(9), 1693-1704 (2004).
3. Pakos-Zebrucka, K., Koryga, I., Mnich, K., *et al.* The integrated stress response. *EMBO Rep.* **17**(10), 1374-1395 (2016).
4. Humeau, J., Leduc, M., Cerrato, G., *et al.* Phosphorylation of eukaryotic initiation factor-2 $\alpha$  (eIF2 $\alpha$ ) in autophagy. *Cell Death Dis.* **11**(6), 433 (2020).
5. Muaddi, H., Majumder, M., Peidis, P., *et al.* Phosphorylation of eIF2 $\alpha$  at serine 51 is an important determinant of cell survival and adaptation to glucose deficiency. *Mol. Biol. Cell* **21**(18), 3220-3231 (2010).