

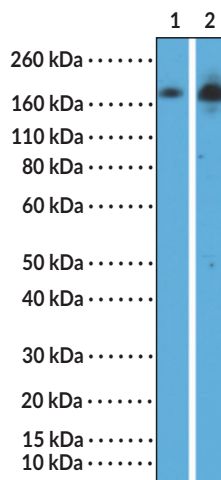
PRODUCT INFORMATION

EGFR (Phospho-Tyr¹⁰⁶⁸) Rabbit Monoclonal Antibody (Clone RM443)
Item No. 35901

Overview and Properties

Contents: This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms: Epidermal Growth Factor Receptor, ErbB-1, HER1
Immunogen: A phospho-peptide corresponding to human phospho-EGFR
Cross Reactivity: (+) EGFR (Phospho-Tyr¹⁰⁶⁸); (-) EGFR without phosphorylation at Tyr¹⁰⁶⁸
Species Reactivity: (+) Human
Form: Liquid
Storage: -20°C (as supplied)
Stability: ≥1 year
Storage Buffer: PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone: RM443
Host: Rabbit
Isotype: IgG
Applications: Western blot (WB); the recommended starting dilution is 1:1,000 -1:4,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: A431 cells nontreated
Lane 2: A431 cells treated

WB of A431 cells nontreated and treated with EGF using EGFR (Phospho-Tyr¹⁰⁶⁸) Rabbit Monoclonal Antibody at a dilution of 1:1,000.

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, 12/07/2021

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

Epidermal growth factor receptor (EGFR), also known as HER1 and ErbB-1, is a cell surface receptor and member of the EGF family of receptor tyrosine kinases with roles in cell proliferation, differentiation, and survival.^{1,2} It is a transmembrane receptor composed of an intracellular tyrosine kinase domain, a transmembrane lipophilic segment, and an extracellular domain that is expressed in epithelial, mesenchymal, and neuronal tissues.¹⁻³ Under unstimulated conditions, EGFR is an auto-inhibited monomer in the plasma membrane.¹ Upon canonical ligand binding, EGFR undergoes homodimerization or heterodimerization with HER2, HER3, or HER4, which induces a conformational change in the cytoplasmic domain that facilitates autophosphorylation and intracellular signaling. EGFR contains five C-terminal autophosphorylation sites, namely tyrosine 1068 (Tyr¹⁰⁶⁸), Tyr¹¹⁴⁸, Tyr¹¹⁷³, Tyr¹⁰⁸⁶, and Tyr⁹⁹².⁴ EGFR (phospho-Tyr¹⁰⁶⁸) is associated with prolonged progression-free survival in patients with non-small cell lung cancer (NSCLC). Cayman's EGFR (Phospho-Tyr¹⁰⁶⁸) Rabbit Monoclonal Antibody can be used for Western blot (WB).

References

1. Sigismund, S., Avanzato, D., and Lanzetti, L. Emerging functions of the EGFR in cancer. *Mol. Oncol.* **12(1)**, 3-20 (2018).
2. Herbst, R.S. Review of epidermal growth factor receptor biology. *Int. J. Radiat. Oncol. Biol. Phys.* **59(2 Suppl)**, 21-26 (2004).
3. Yano, S., Kondo, K., Yamaguchi, M., *et al.* Distribution and function of EGFR in human tissue and the effect of EGFR tyrosine kinase inhibition. *Anticancer Res.* **23(5A)**, 3639-3650 (2003).
4. Wang, F., Wang, S., Wang, Z., *et al.* Phosphorylated EGFR expression may predict outcome of EGFR-TKIs therapy for the advanced NSCLC patients with wild-type EGFR. *J. Exp. Clin. Cancer Res.* **31(1)**, 65 (2012).

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