

# PRODUCT INFORMATION

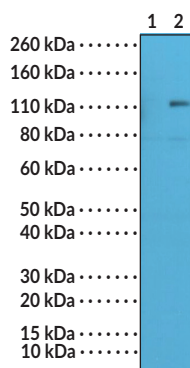


## EGFR (Phospho-Tyr<sup>1173</sup>) Rabbit Monoclonal Antibody (Clone RM269) Item No. 32219

### Overview and Properties

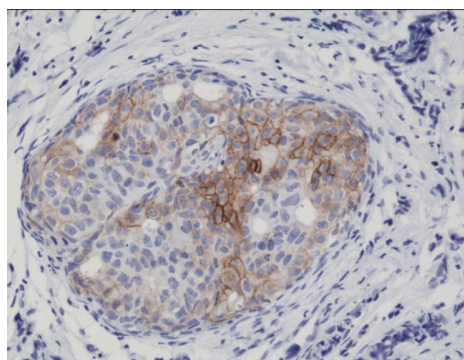
<b>Contents:</b>	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	Epidermal Growth Factor Receptor, ErbB-1, HER1
<b>Immunogen:</b>	Peptide corresponding to human EGFR (phospho-Tyr <sup>1173</sup> )
<b>Cross Reactivity:</b>	(+) EGFR (phospho-Tyr <sup>1173</sup> ); (-) EGFR without phosphorylation at Tyr <sup>1173</sup>
<b>Species Reactivity:</b>	(+) Human
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM269
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:500-1:1,000 and 1:1,000-1:5,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: A431 cells (untreated)  
Lane 2: A431 cells (treated)

WB of A431 cells untreated or treated with EGF using EGFR (Phospho-Tyr<sup>1173</sup>) Rabbit Monoclonal Antibody (Clone RM269) at a dilution of 1:500.



Immunohistochemical staining of formalin-fixed, paraffin-embedded (FFPE) human breast cancer tissue sections using EGFR (Phospho-Tyr<sup>1173</sup>) Rabbit Monoclonal Antibody at a dilution of 1:5,000.

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, 02/05/2024

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

---

Epidermal growth factor receptor (EGFR), also known as HER1 and ERBB1, is a cell surface receptor and member of the EGF family of receptor tyrosine kinases with roles in cell proliferation, differentiation, and survival.<sup>1,2</sup> 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.<sup>1-3</sup> Under unstimulated conditions, EGFR is an auto-inhibited monomer in the plasma membrane.<sup>1</sup> 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, tyrosine 1068 (Tyr<sup>1068</sup>), Tyr<sup>1148</sup>, Tyr<sup>1173</sup>, Tyr<sup>1086</sup>, and Tyr<sup>992</sup>.<sup>4</sup> EGFR autophosphorylation at Tyr<sup>1173</sup> (phospho-Tyr<sup>1173</sup>) leads to interaction with phospholipase C $\gamma$  (PLC $\gamma$ ) and Shc, which have roles in activating the MAPK signaling pathway. Increased levels of EGFR (phospho-Tyr<sup>1173</sup>) are associated with poor progression-free survival in patients with non-small cell lung cancer (NSCLC). Levels of EGFR (phospho-Tyr<sup>1173</sup>) are increased in, and positively correlate with activation of MAPK signaling and cytokine production in, bronchial epithelial biopsies from healthy individuals exposed to diesel exhaust.<sup>5</sup> Cayman's EGFR (Phospho-Tyr<sup>1173</sup>) Rabbit Monoclonal Antibody (Clone RM269) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

## 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).
5. Pourazar, J., Blomberg, A., Kelly, F.J., et al. Diesel exhaust increases EGFR and phosphorylated C-terminal Tyr 1173 in the bronchial epithelium. *Part. Fibre Toxicol.* **5**, 8 (2008).

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