

PRODUCT INFORMATION

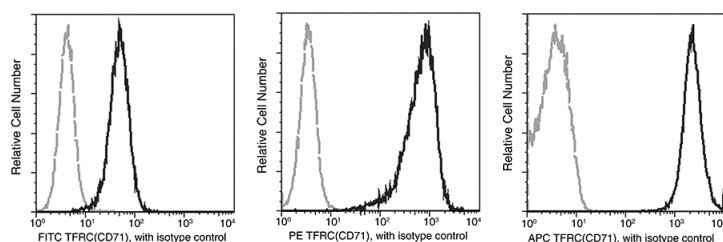


Transferrin Receptor Protein 1/CD71 Monoclonal Antibody (PE) (Clone 04) Item No. 37081

Overview and Properties

| | |
|----------------------------|--|
| Contents: | 250 µl or 1 ml of protein A-affinity purified monoclonal antibody. |
| Synonyms: | Cluster of Differentiation 71, p90, TfR1, TFRC |
| Immunogen: | Recombinant human TfR1 |
| Cross Reactivity: | (+) TfR1 |
| Species Reactivity: | (+) Human |
| Form: | Liquid |
| Storage: | 2-8°C (as supplied) |
| Stability: | ≥1 year |
| Storage Buffer: | PBS with 0.5% BSA and 0.09% sodium azide |
| Concentration: | 100 µg/ml |
| Clone: | 04 |
| Host: | Mouse |
| Isotype: | IgG1 |
| Application: | Flow Cytometry; the optimal working concentration/dilution should be determined empirically. |

Images



Flow cytometric analysis of human TfR1/CD71 expression on U937 cells. Cells were stained with Transferrin Receptor Protein 1/CD71 Monoclonal Antibody (PE) (Clone 04). The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

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, 02/07/2024

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

Transferrin receptor (TfR1), also known as CD71, is a homodimeric transmembrane receptor for transferrin (Item No. 32030) that facilitates iron delivery into cells and is encoded by *TFRC* in humans.¹ It is composed of two TfR1 monomers, each containing a cytoplasmic tail, an internalization motif, a membrane-spanning portion, and a stalk region that covalently links the monomers. An extracellular ectodomain binds transferrin and drives TfR1 dimerization.^{1,2} TfR1 is ubiquitously expressed, except on mature red blood cells and certain terminally differentiated cells, with the highest expression on immature erythroid cells and in the placenta, and is involved in erythropoiesis, lymphocyte development, and hematopoietic expansion in the bone marrow.^{3,4} TfR1/transferrin-mediated iron transport contributes to the intracellular iron pool required for ferroptosis and an anti-TfR1 antibody has been used in combination with anti-malondialdehyde antibodies to identify ferroptotic cells *in vitro* and human cancer tissue in a mouse xenograft model.⁵ *TFRC* is overexpressed in various breast cancer tumors and gliomas and positively correlated with poor prognosis.⁶ Cayman's Transferrin Receptor Protein 1/CD71 Monoclonal Antibody (PE) (Clone 04) is composed of a TfR1 monoclonal antibody conjugated to phycoerythrin (PE) and can be used for flow cytometry.

References

1. Luck, A.N. and Mason, A.B. Transferrin-mediated cellular iron delivery. *Curr. Top. Membr.* **69**, 3-35 (2012).
2. Gammella, E., Buratti, P., Cairo, G., *et al.* The transferrin receptor: The cellular iron gate. *Metallomics* **9**(10), 1367-1375 (2017).
3. Ponka, P. and Lok, C.N. The transferrin receptor: Role in health and disease. *Int. J. Biochem. Cell Biol.* **31**(10), 1111-1137 (1999).
4. Wang, S., He, X., Wu, Q., *et al.* Transferrin receptor 1-mediated iron uptake plays an essential role in hematopoiesis. *Haematologica* **105**(8), 2071-2082 (2020).
5. Feng, H., Schorpp, K., Jin, J., *et al.* Transferrin receptor is a specific ferroptosis marker. *Cell Rep.* **30**(10), 3411-3423 (2020).
6. Shen, Y., Li, X., Dong, D., *et al.* Transferrin receptor 1 in cancer: A new sight for cancer therapy. *Am. J. Cancer Res.* **8**(6), 916-931 (2018).

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