

CFTR monoclonal antibody (CF3)

Cystic Fibrosis (CF) is a common lethal genetic disease caused by mutations of the gene coding for the cystic fibrosis transmembrane conductance factor, a cAMP regulated chloride channel. Approximately 70% of all CF cases share the deletion of a phenylalanine at position 508 (δ F508) which results in abnormal chloride transport. Since the CF mutation is lethal, most often by lung and liver disease, it raises the question of why this genetic disease remains as common as it is. One possible explanation is that *Salmonella typhi* has been shown to use CFTR to enter intestinal epithelial cells and that delta F508 heterozygote and homozygote mice showed 86% and 100% reductions in *S. typhi* intestinal submucosal uptake.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cystic fibrosis transmembrane conductance regulator, ABCC7
Application	ICC, IP, WB
Application Notes	Detects a band of ~170kDa by Western blot. Functional Application: inhibits epithelial uptake of <i>S. typhi</i> in some mouse cell lines.
Clone	CF3
Crossreactivity	Cross-reacts with related proteins in mouse cell line Heb7a that doesn not contain CFTR mRNA.
Formulation	Liquid. Diluted ascites containing 0.05% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 103-117(G103RIIASYDPDNKEER117) in the first extracellular loop of human and rabbit CFTR (cystic fibrosis transmembrane conductance regulator). This sequence is highly conserved in mouse, sheep, bovine and <i>Xenopus laevis</i> CFTR.
Isotype	IgM

Recommendation Dilutions/Conditions

Immunocytochemistry (1:500; staining of CFTR in mouse epithelial cells results in cell surface staining, consistent with localization at the plasma membrane. Antibody also detects one or more immunologically related proteins in murine cell line Heb7a that does not contain CFTR mRNA.)Western Blot (1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity

Human, Mouse

UniProt ID

P13569

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com