

Cytokeratin 7, 17 & 19 monoclonal antibody (4F5)

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

Ordering Information

[Order Online »](#)

BML-CA1139-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CK7/CK17/CK19
Application	ICC, WB
Application Notes	Detects bands of 54 kDa, 46 kDa and 40 kDa by Western Blot.
Clone	4F5
Formulation	Liquid. Containing 0.1% sodium azide.
Host	Mouse
Immunogen	Cytokeratins isolated from cultured porcine kidney epithelial cells (LLC-PK).
Isotype	IgG1
Purity Detail	Protein A affinity chromatography.
Species Reactivity	Human, Mouse, Porcine, Rat
Specificity	Recognizes cytokeratin proteins 7, 17 and 19.
UniProt ID	P08729 (human CK7), Q04695 (human CK17), P08727 (human CK19)

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
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@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