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 NotesDetects 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).

lgG1

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

