Cytokeratin 18 monoclonal antibody (4B11)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-CA1138-0100

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 CK18

Application ICC, WB

Clone 4B11

Host Mouse

Immunogen Cytokeratins isolated from cultured porcine kidney

epithelial cells (LLC-PK).

lsotype lgG1

Purity Detail Purified.

Species Reactivity Human, Mouse, Porcine, Rat

Specificity Recognizes cytokeratin 18 in a variety of species including

human, mouse, rat and pig.

UniProt ID P05783 (human)

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

