# Plk1 monoclonal antibody (3F8)

The polo-like kinases (PLKs) function in cell-cycle progression and in multiple stages of mitosis. PLK1 is activated by phosphorylation of Thr210 at the G2/M phase boundary. Active PLK1 participates in mitotic entry through activation of Cdc25C and nuclear import of cyclin B1, thereby activating the Cdc2/cyclin B kinase.

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

**View Online »** 

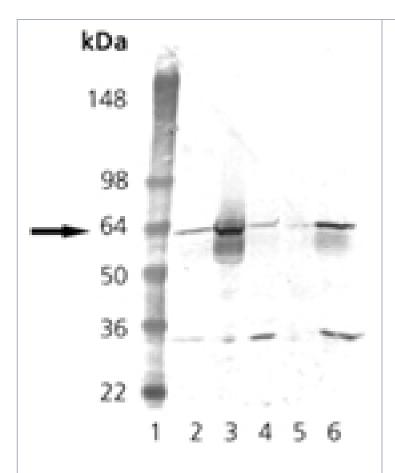
## **Ordering Information**

**Order Online »** 

ADI-KAM-CC105-D	50µg
ADI-KAM-CC105-F	200μg

Manuals, SDS & CofA

**View Online »** 



Western blot analysis of PLK1 mAb (3F8): Lane 1: MW marker, Lane 2: HeLa, Lane 3: Jurkat, Lane 4: PC-12, Lane 5: 3T3, Lane 6: HEK-293.

### **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 Polo-like kinase 1

**Application** WB

**Application Notes**Detects a band of ~62kDa by Western blot.

Clone 3F8

**Crossreactivity** Weakly cross-reacts with mouse Plk1.

**Formulation** Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09%

sodium azide.

**Host** Mouse

**Immunogen** Recombinant fragment from the middle of human PLK1.

lsotype IgG1

Purity Detail Protein G affinity purified.

**Recommendation Dilutions/Conditions**Western Blot (1:1,000, colorimetric)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

**Source** Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Rat

UniProt ID P53350

Last modified: May 29, 2024



info-

eu@enzolifesciences.com