

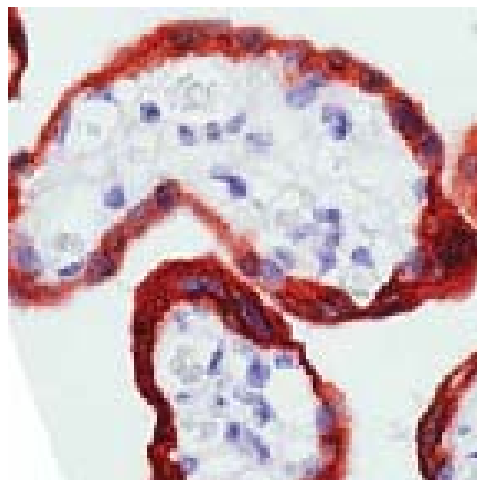


## Product Data Sheet

Product Name:	Anti - c - Abl (monoclonal)
Catalog Number:	53278
Lot Number:	See label on vial
Product Description:	This mouse monoclonal antibody, purified from ascites fluid by protein G, is supplied as 50 µg in 250 µl (200 µg/ml) of 1X PBS containing with 0.2% BSA and 0.05% sodium azide.
Immunogen:	Mouse Anti-c-ABL monoclonal antibody was raised against a recombinant ABL protein with Ig isotype being IgG <sub>1</sub> /k.
Species Reactivity:	Species reactivity includes human and mouse, while others remain unknown.
Application Notes:	The following concentration ranges are recommended starting points for this product. Optimal working concentrations should be determined by the investigator for specific applications.

WB:	2.5-5.0 µg/ml
IF:	Not Tested
IP:	2.0 µg/mg protein lysate
IHC:	1:50 for 30 min at room temperature

Staining of formalin-fixed tissues requires boiling tissue sections in 1 mM EDTA, pH 8.0, for 10 min followed by cooling at room temperature for 20 min.



**Left:** Human placenta stained with anti-c-ABL

**Positive Control:**  
HeLa cells or placenta

**Cellular Localization:**  
Cytoplasmic and nuclear

**Background:**

The c-ABL proto-oncogene encodes a protein tyrosine kinase that is located in the cytoplasm and nucleus. In chronic myelogenous leukemia and in a subset of acute lymphoblastic leukemias, the c-ABL proto-oncogene undergoes a chromosomal translocation producing a novel rearranged chromosome (the Philadelphia chromosome) as a result of the fusion of c-ABL sequences from chromosome 9 to the Bcr gene on chromosome 22. The molecular consequence of this translocation is the generation of a chimeric Bcr/ABL mRNA encoding activated ABL protein tyrosine kinase.

**Storage:**

Store at 2-8 °C for up to one year. Avoid repeated freezing and thawing.

This product is for *in vitro* research use only.