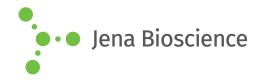
DATA SHEET

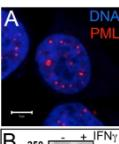


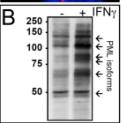


anti-PML anti-human PML rabbit, polyclonal

Cat. No. **Amount**

ABD-030 50 μl





For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles, store in aliquots

Shelf Life: 12 months

Form: liquid (Supplied as serum)

Applications:

Indirect Immunofluorescence 1:250. Western Blot 1:500

Description:

Polyclonal Antibody pABahPML(460-479)#3 to human PML.

Antibody specificity was evalutated by indirect immunoflourescence and western blotting (see Figure 1A and B). The antibody detects all PML isoforms according to Condemine et al. (2006).

Immunogen:

The antibody binds to a peptide corresponding to amino acids 460-479 (KGPSYGEDVSNTTTQKRKC) of the human PML protein. No cross reaction with mouse PML protein was observed in Western blots. The immunogen peptide sequence is not conserved in rat, chimpanzee, dog, or bovine PML proteins.

Selected References:

Krejcí et al. (2008) Nuclear organization of PML bodies in leukaemic and multiple myeloma cells. Leukemia Research 32:1866.

Condemine et al. (2006) Characterization of endogenous human promyelocytic leukemia isoforms. Cancer Res. 66:6192.