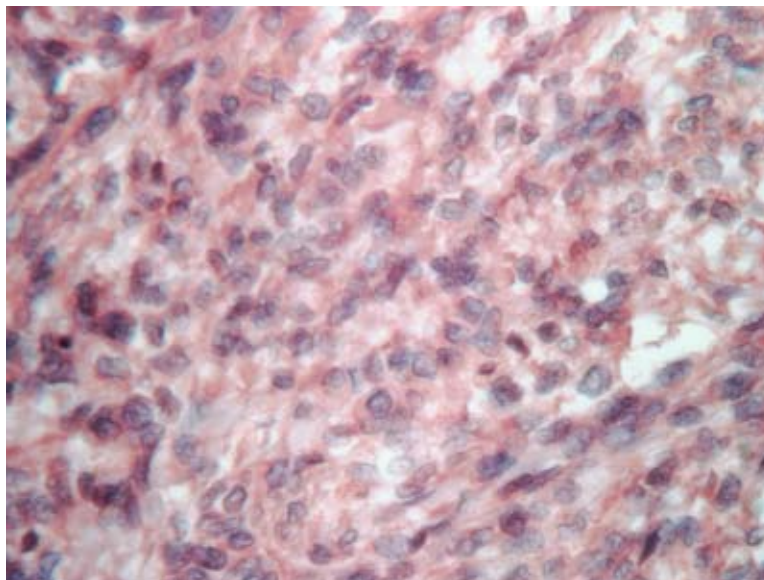




Product Data Sheet

Product Name:	Anti - CD117/c - kit
Catalog Number:	29562
Lot Number:	See label on vial
Product Description:	This polyclonal antibody is supplied as an epitope affinity purified rabbit IgG, 50 µg in 250 µl of phosphate buffered saline (pH 7.4) containing 0.02% sodium azide.
Immunogen:	Rabbit anti-CD117 polyclonal antibody was raised against a synthetic peptide corresponding to tail portion of human CD117/c-kit.
Species Reactivity:	This antibody is specific to human and dog.
Application Notes:	The following concentration ranges are recommended starting points for this product. Western blot: 0.5 - 2 µg/ml IHC: 1:400 for 10 min @ RT



Above: Human Stroma tumor stained with anti-CD117 (Cat# 29562) by IHC Histain Kit-AEC (Cat# 29702).

Background:

The CD117 molecule, a tyrosine kinase receptor with a molecular weight of 145 kDa, is encoded by proto-oncogene c-kit (1). The stem cell factor (SCF) binding to the CD117/c-kit receptor triggers homodimerization and intermolecular tyrosine phosphorylation, thus initiating signal transduction (2). The receptor tyrosine kinase CD117/c-kit and its ligand SCF are essential to the maturation of hemopoietic, primordial germ cells precursors and melanocytes during embryonic development. Multiple cellular functions are affected by CD117/c-kit dependent signals including cell survival, proliferation, adhesion, differentiation and functional maturation of hemopoietic cells (3-4).

Storage:

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

References:

1. Elmore, L. et al. *Arch Pathol Lab Med* **125**, 146 (2001).
2. Ahman, L. et al. *Leuk Lymphoma* **34**, 451 (1999).
3. Broudy, V. et al. *Blood* **91**, 898 (1998).
4. Alexander, W. et al. *Leuk Res* **22**, 1097 (1998).

This product is for in vitro research use only.