



EAGLE
BIOSCIENCES

Anti-Rabbit Cytochrome b5

Product Number: CB500-A100

Aliquot: 100 μ L

Storage: -80°C

PRODUCT SPECIFICATION SHEET

DESCRIPTION:	This is a goat polyclonal antibody and is specific for cytochrome b5 in rabbit but cross reacts with human and rat. This antibody is non-cross reactive with the 50 and 60 kDa species of cytochrome b5.
CONCENTRATION:	Not determined.
PRESENTATION:	Neat antiserum with 0.08% (w/v) sodium azide.
STORAGE:	-80°C. AVOID REPEATED FREEZE THAWS.
IMMUNOGEN:	Rabbit cytochrome b5
PURIFICATION:	Not purified
APPLICATIONS:	This antibody is suitable for use in Western blotting at 1:1000 dilution.
SPECIES SPECIFICITY:	Rabbit, Rat and Human.

Product Developed and Manufactured in the USA

EAGLE BIOSCIENCES, INC.
20A NW Blvd., Suite 112
Nashua, NH 03063

Phone: 617-419-2019
Fax: 617-419-1110
Email: info@EagleBio.com
www.EagleBio.com