

## SCIENTIFIC BACKGROUND:

Choline acetyltransferase is found in cholinergic neurons on brain and central nervous system.

## SPECIFICITY:

This anti-choline acetyltransferase antibody recognizes choline acetyltransferase found in cholinergic neurons in the brain and central nervous system of rat, mouse, human, guinea pig, chicken, opossum, avian and macaque monkey (immunoblot).

## IMMUNOGEN:

This goat polyclonal antibody was generated using human placental choline acetyltransferase enzyme.

## FORMAT:

0.5 mL affinity purified goat polyclonal antibody in PBS containing 5 mg/mL BSA with 0.2% sodium azide.

## STABILITY AND STORAGE:

Store at -20°C in undiluted aliquots. Under these conditions, product is stable for up to 12 months from date of receipt.

**Avoid repeated freezing and thawing.**

## APPLICATIONS AND DIRECTIONS FOR USE:

Centrifuge tube briefly before use to ensure recovery of entire content.

Dilute with medium or phosphate buffered saline containing appropriate blocking serum. The suggested working dilution for immunoblot is 1:100-1,000. This antibody reacts with a band approximately 68-70 kDa. For immunohistochemistry, the suggested working dilution is 1:100. The optimal working dilution should be determined for each specific application.

**THIS REAGENT IS FOR RESEARCH ONLY.  
IT IS NOT TO BE ADMINISTERED TO HUMANS.**

**Hazardous Ingredient: Sodium Azide.** Avoid exposure to skin and eyes, ingestion and contact with heat, acids and metals. Wash exposed skin with soap and water. Flush eyes with water. Dilute with running water before discharging into plumbing.


Product Information Sheet

**ANTI-CHOLINE  
ACETYLTRANSFERASE  
(ChAT)**

Affinity Purified Goat Polyclonal  
Antibody

Catalog # 01510 0.5 mL

**ANTIBODIES**



## StemCell Technologies

### In North America

Tel: 1.604.877.0713  
Fax: 1.604.877.0704  
Toll Free Tel: 1.800.667.0322  
Toll Free Fax: 1.800.567.2899  
e-mail: [info@stemcell.com](mailto:info@stemcell.com)  
[www.stemcell.com](http://www.stemcell.com)

### In the United Kingdom

Tel: +44.(0).20.7537.7565  
Fax: +44.(0).20.7515.5408  
Toll Free within United Kingdom  
Tel: 0800.731.27.14  
Fax: 0800.731.27.13  
e-mail: [info@stemcellgb.com](mailto:info@stemcellgb.com)  
[www.stemcellgb.com](http://www.stemcellgb.com)

### In Europe

Tel: +33.(0).4.76.04.75.30  
Fax: +33.(0).4.76.18.99.63  
e-mail: [info@stemcellfrance.com](mailto:info@stemcellfrance.com)

Revised:  
October 2004