

## SCIENTIFIC BACKGROUND:

The anti-GFAP antibody can be used to detect glial fibrillary acidic protein, a differentiation marker for astrocytes, localized in normal central nervous system tissue and certain tumor and metastases of the glial antigen. The antibody is also suitable for immunofluorescent labeling of cultured mammalian cells.

## SPECIFICITY:

This antibody reacts with glial fibrillary acidic protein (GFAP).

## ISOTYPE:

IgG fraction of rabbit antiserum.

## IMMUNOGEN:

The antiserum was developed in rabbits using GFAP purified from human brain as the immunogen. The whole antiserum was fractionated and purified by ion-exchange chromatography to isolate the IgG fraction which is essentially free of rabbit serum proteins.

## FORMAT:

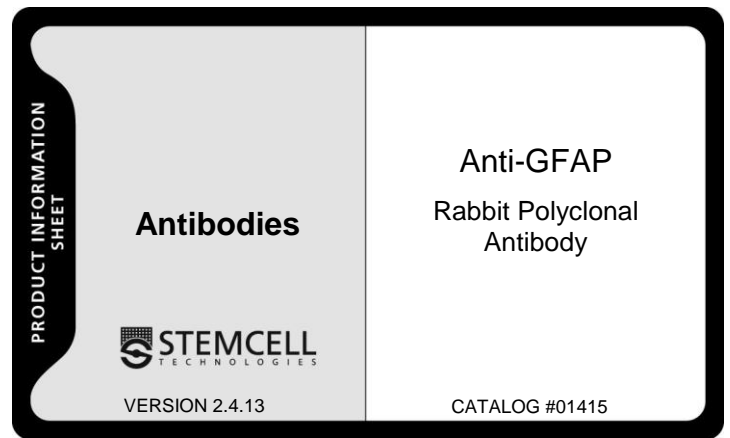
200 µL in 0.01 M phosphate buffered saline, pH 7.4, with 15 mM sodium azide.

**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.

**Concentration will vary from lot to lot.** Please refer to Table 1 for lot-specific concentration. If your lot number differs from those shown, please contact STEMCELL Technologies for information specific to your antibody.

Table 1. Lot-Specific Antibody Concentrations

LOT#	CONCENTRATION
11K41730	9.1 mg/mL
11G40408	9.1 mg/mL
11F40089	9.1 mg/mL
11F39909	9.1 mg/mL
11B38532	9.1 mg/mL
11A38037	9.1 mg/mL
10M37819	9.1 mg/mL
10J36413	9.1 mg/mL
10G35698	9.1 mg/mL
10B33705	9.1 mg/mL
10A33095	9.1 mg/mL
09L32286	9.1 mg/mL



## STABILITY AND STORAGE:

Store at 2 - 8°C in undiluted aliquots for up to one month. For extended storage, store at -20°C in undiluted aliquots. Repeated freezing and thawing is not recommended.

## APPLICATIONS AND DIRECTIONS FOR USE:

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

Dilute with medium of phosphate buffered saline containing appropriate blocking serum. The optimal working dilution should be determined for each specific application. Refer to the NeuroCult® Procedure Manual or ES-Cult® Protocols – Neural available on our website at [www.stemcell.com](http://www.stemcell.com) for further instructions on how to use this antibody.