

# PRODUCT INFORMATION



## GFAP (human, recombinant)

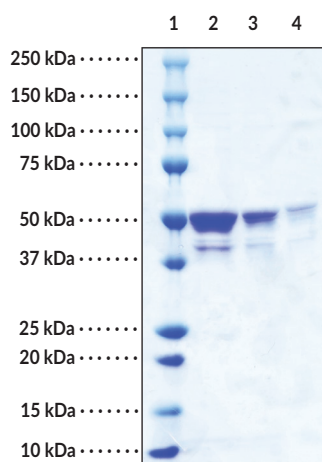
Item No. 27353

### Overview and Properties

**Synonyms:** ALXDRD, Glial Fibrillary Acidic Protein, Intermediate Filament Protein  
**Source:** Recombinant N-terminal histidine-tagged GFAP expressed in *E. coli*  
**Amino Acids:** 2-432  
**Uniprot No.:** P14136  
**Molecular Weight:** 51.9 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** *batch specific* (≥65% estimated by SDS-PAGE)  
**Supplied in:** 10 mM Tris, pH 8.0  
**Protein**  
**Concentration:** *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: GFAP (8 µg)  
Lane 3: GFAP (4 µg)  
Lane 4: GFAP (2 µg)

Representative gel image shown; actual purity may vary between each batch.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/17/2020

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

Glial fibrillary acidic protein (GFAP) is a protein encoded by the *GFAP* gene in humans and a member of the class III intermediate filament (IF) protein family.<sup>1</sup> It is composed of an N-terminal head domain, a highly conserved  $\alpha$ -helical rod domain, and a C-terminal tail domain that mediate GFAP self-assembly, dimerization, and oligomerization, respectively.<sup>2,3</sup> GFAP is expressed in, and has commonly been used as a pan marker for, mature astrocytes.<sup>1</sup> GFAP IFs form a dynamic network of cytosolic filament proteins that collectively provide structure and strength to the cytoskeleton of astrocytes, thus supporting their morphology and function.<sup>1</sup> Isolated astrocytes from neonatal *Gfap*<sup>-/-</sup> mouse brain have reduced numbers of IFs and IF bundles, increased proliferation, and loss of contact-inhibited growth.<sup>4,5</sup> *Gfap*<sup>-/-</sup> mice develop more diffuse and infiltrative brain lesions compared to wild-type littermates in a mouse model of experimental autoimmune encephalomyelitis (EAE).<sup>6</sup> Mutations in the rod and tail domains of *GFAP* have been associated with Rosenthal fiber formation, a hallmark of Alexander disease.<sup>7</sup> Transgenic overexpression of *Gfap* in mice increases the expression of certain cytokines and antioxidative enzymes in the olfactory bulb and has been used as a mouse model of Alexander disease.<sup>8</sup> GFAP can be citrullinated on the arginine residue at position 270 (R270) and at R416 by protein arginine deiminase 1 (PAD1; Item No. 10784) and PAD2 (Item No. 10785).<sup>9</sup> Citrullinated GFAP has been found in rat cerebral cortex in a model of traumatic brain injury, as well as in postmortem hippocampus from patients with Alzheimer's disease.<sup>9,10</sup>

## References

1. Hol, E.M. and Capetanaki, Y. Type III intermediate filaments desmin, glial fibrillary acidic protein (GFAP), vimentin, and peripherin. *Cold Spring Harb. Perspect. Biol.* **9**(12), a021642 (2017).
2. Inagaki, M., Nakamura, Y., Takeda, M., et al. Glial fibrillary acidic protein: Dynamic property and regulation by phosphorylation. *Brain Pathol.* **4**(3), 239-243 (1994).
3. Chen, W.-J. and Liem, R.K.H. The endless story of the glial fibrillary acidic protein. *J. Cell Sci.* **107**(Pt 8), 2299-2311 (1994).
4. Pekny, M., Eliasson, C., Chien, C.L., et al. GFAP-deficient astrocytes are capable of stellation *in vitro* when cocultured with neurons and exhibit a reduced amount of intermediate filaments and an increased cell saturation density. *Exp. Cell Res.* **239**(2), 332-343 (1998).
5. Rutka, J.T. and Smith, S.L. Transfection of human astrocytoma cells with glial fibrillary acidic protein complementary DNA: Analysis of expression, proliferation, and tumorigenicity. *Cancer Res.* **53**(15), 3624-3631 (1993).
6. Liedtke, W., Edelmann, W., Chiu, F.C., et al. Experimental autoimmune encephalomyelitis in mice lacking glial fibrillary acidic protein is characterized by a more severe clinical course and an infiltrative central nervous system lesion. *Am. J. Pathol.* **152**(1), 251-259 (1998).
7. Li, R., Messing, A., Goldman, J.E., et al. GFAP mutations in Alexander disease. *Int. J. Dev. Neurosci.* **20**(3-5), 259-268 (2002).
8. Hagemann, T.L., Gaeta, S.A., Smith, M.A., et al. Gene expression analysis in mice with elevated glial fibrillary acidic protein and Rosenthal fibers reveals a stress response followed by glial activation and neuronal dysfunction. *Hum. Mol. Genet.* **14**(16), 2443-2458 (2005).
9. Ishigami, A., Masutomi, H., Handa, S., et al. Mass spectrometric identification of citrullination sites and immunohistochemical detection of citrullinated glial fibrillary acidic protein in Alzheimer's disease brains. *J. Neurosci. Res.* **93**(11), 1664-1674 (2015).
10. Lazarus, R.C., Buonora, J.E., Flora, M.N., et al. Protein citrullination: A proposed mechanism for pathology in traumatic brain injury. *Front. Neurol.* **6**:204 (2015).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM