

S100A1

Calcium-binding Protein A1

human, recombinant, *E. coli*

Cat. No.	Amount
PR-391	50 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied as 100 μ g/ml solution in 20 mM HEPES-KOH pH 7.0, 50 mM NaCl, 1 mM EDTA and 1 mM DTT.

Activity

Assayed for calcium binding.

Purity

>95% by SDS-PAGE.

Description

The proteins of the S100 family belong to the family of calcium binding proteins with EF-hand type Ca^{2+} binding motive. S100A1 (calcium-binding protein A1) is involved in the activation of sarcoplasmic calcium release and the regulation of intermediate filament polymerisation.

Selected References:

Zimmer *et al.* (1995) The S100 protein family - history, function, and expression. *Brain Research Bulletin* **37**:417.

Heizmann *et al.* (1998) New perspectives on S100 proteins: a multi-functional Ca^{2+} , Zn^{2+} , and Cu^{2+} -binding protein family. *Biometals* **11**:383.

Arcuri *et al.* (2002) Annexin V, Annexin VI, S100A1 and S100B in developing and adult avian skeletal muscles. *Neuroscience* **109**:371.