

## S100A1

**Calcium-binding Protein A1**  
human, recombinant, *E. coli*

Cat. No.	Amount
PR-391	50 µg

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

### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 20 mM Tris-HCl pH 7.0,  
0.1 mM EDTA.

#### 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<sup>2+</sup> 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<sup>2+</sup>-, Zn<sup>2+</sup>-, and Cu<sup>2+</sup>-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.