

S100A11

Calcium-binding Protein A11

human, recombinant, *E. coli*

Cat. No.	Amount
PR-1265	10 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Protein synonyms/aliases

Calgizzarin (S100 calcium-binding protein A11),
S100C protein and MLN70.

Form

Liquid. Supplied in PBS and 50% glycerol.

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 Ca²⁺ binding motive.

S100A is composed of an alpha and beta chain whereas S100B is composed of two beta chains. S100A11 is implicated in a variety of biologic functions such as proliferation and differentiation as well as in cancer. S100A11 expression is associated with the development of lymph node metastases and appears to be a molecular marker for clinical prognosis.

Recombinant Human S100A11 is expressed in *E. coli* having a molecular weight of 17 kDa fused to an amino terminal His tag.

Selected References:

- Pan *et al.* (2006) Folding Kinetics of the S100A11 Protein Dimer Studied by Time-Resolved Electrospray Mass Spectrometry and Pulsed Hydrogen-Deuterium Exchange. *Biochemistry*. **45**:3005.
- Rust *et al.* (2005) High expression of calcium-binding proteins, S100A10, S100A11 and CALM2 in anaplastic large cell lymphoma. *Br J Haematol*. **131**:596.
- Melle *et al.* (2006) Different expression of calgizzarin (S100A11) in normal colonic epithelium, adenoma and colorectal carcinoma. *Int J Oncol*. **28**:195.
- Cross *et al.* (2005) Expression of S100 proteins in normal human tissues and common cancers using tissue microarrays: S100A6, S100A8, S100A9 and S100A11 are all overexpressed in common cancers. *Histopathology*. **46**:256.
- Rehman *et al.* (2004) Dysregulated expression of S100A11 (calgizzarin) in prostate cancer and precursor lesions. *Hum Pathol*. **35**:1385.
- Mori *et al.* (2004) S100A11 gene identified by in-house cDNA microarray as an accurate predictor of lymph node metastases of gastric cancer. *Oncol Rep*. **11**:1287.