

## S100A6

### Calcium-binding Protein A6

human, recombinant, *E. coli*

Cat. No.	Amount
PR-1260	5 µg

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

#### Avoid freeze / thaw cycles

#### Protein synonyms/aliases

S100 calcium-binding protein A6 and Calcyclin.

#### Form

Lyophilized (no additives).

#### Solubility

It is recommended to reconstitute the lyophilized PDGF-AB in sterile bidest H<sub>2</sub>O not less than 100 µg/ml. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).

#### Activity

Assayed for calcium binding.

#### Application

May be used in Immunoblots, Immunohistochemistry and for Absorption.

#### Purity

>90% by SDS-PAGE (reducing and non reducing)

#### Description

The proteins of the S100 family belongs to the family of calcium binding proteins with EF-hand type Ca<sup>2+</sup> binding motive. S100A is composed of an alpha and beta chain whereas S100B is composed of two beta chains.

S100A6 (calcium-binding protein A6) is involved in the invasive process of human carcinomas (e.g. colorectal adenocarcinomas, pancreatic and prostate cancer). S100A6 expression levels decrease when carcinoma cells form glandular structures at the central portions of metastatic nodules.

S100A6 appears to be a marker in clinical cancer diagnostic.

The protein is purified by HPLC.

#### Selected References:

Filipek (2006) S100A6 and CacyBP/SIP - two proteins discovered in ehrlich ascites tumor cells that are potentially involved in the degradation of beta-catenin. *Chemotherapy*. **52**:32.

Cheng *et al.* (2005) Calcium-binding proteins annexin A2 and S100A6 are sensors of tubular injury and recovery in acute renal failure. *Kidney Int.* **68**:2694.

Lesniak *et al.* (2005) Binding and functional characteristics of two E-box motifs within the S100A6 (calcyclin) gene promoter. *J Cell Biochem.* **97**(5):1017.

Ohuchida *et al.* (2005) The role of S100A6 in pancreatic cancer development and its clinical implication as a diagnostic marker and therapeutic target. *Clin Cancer Res.* **11**:7785.

Luu *et al.* (2005) Increased expression of S100A6 is associated with decreased metastasis and inhibition of cell migration and anchorage independent growth in human osteosarcoma. *Cancer Lett.* **229**:135.

Rehman *et al.* (2005) Promoter hyper-methylation of calcium binding proteins S100A6 and S100A2 in human prostate cancer. *Prostate.* **65**:322.