
Anti-p66 α

(Catalog # A-1015)

Background:	Silencing by methylated DNA has been shown to occur when methyl-CpG-binding proteins (MeCP) recruit histone deacetylase complexes. The MeCP1 (methyl-CpG-binding protein 1) complex includes the zinc finger-containing protein known as p66. The p66 α polyclonal antibody detects p66 α , which is thought to play a role in transcription repression.
Concentration:	0.5 mg/ml
Description:	Rabbit polyclonal antibody to p66 α
Purification:	Protein A Purified
Specificity:	Mouse and human
Isotype:	IgG
Formulation:	0.02% Sodium Azide, 0.1 mg/ml BSA, PBS
Storage:	4°C, stable for 6 months from the date of shipment. For long-term storage, aliquot and store at -20°C.
Application:	Western blot: 1:200-1:500 IHC: 1:100 ELISA: 1:1000-1:2000
Research Use:	Research use only.
Handling Recommendations:	For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Ordering Information

Products	Size	Cat. No.
Anti-p66 α	50 μ g	A-1015-050
	100 μ g	A-1015-100

This product is for research purposes only. Not intended for use in diagnostic procedures.