# [pSer<sup>36</sup>]Shc monoclonal antibody (6E10)

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

**Ordering Information** 

Order Online »

ALX-804-358-C100

100µg

Manuals, SDS & CofA

**View Online »** 

### **Handling & Storage**

**Use/Stability** Stable at -80°C up to 1 year, at 4°C up to 3 months.

**Handling** Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid

nitrogen.

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Shc transforming protein 1, SH2 domain protein C1, Shc1

**Application** WB

Clone 6E10

**Crossreactivity** Does not cross-react with non-phosphorylated shc.

Formulation Lyophilized from 1ml of 2x PBS containing 0.09% sodium

azide, PEG, and sucrose.

**Host** Mouse

Immunogen Synthetic peptide corresponding to aa 33-39

(E33LPpSPSA39) of shc phosphorylated at Ser36

conjugated to KLH.

lsotype lgG1

Positive Control Included. (Prod. No. 840-033-1)

Purity Detail Thiophilic adsorption and size exclusion chromatography

purified.

#### **Recommendation Dilutions/Conditions**

Western Blot (1µg/ml for HRPO/ECL detection; recommended blocking buffer BPPT: 1% (w/v) BSA, 1% (w/v) PEG 4,000, 1% (w/v) polyvinylpyrrolidone, 0.1% (v/v) Tween 20, 2 x PBS). Immunoprecipitation of shc prior to detection with MAb to Shc (phosphorylated) (pSer36) (6E10) is recommendedSuggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Reconstitution

Reconstitute with 1ml water (15 minutes at room temperature).

**Species Reactivity** 

Dog, Human, Mouse

**UniProt ID** 

P29353

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024