
SHC Polyclonal Antibody

(Catalog No. A-0494)

Background

There are 3 alternatively spliced isoforms of the Shc adapter protein: p46SHC, p52SHC and p66SHC, that link activated growth factor receptors to signalling pathway. Isoform p46SHC and p52SHC, once phosphorylated, couple activated receptor tyrosine kinases to Ras via the recruitment of the GRB2/SOS complex and are implicated in mitogenic signalling. They function as initiators of the Ras signalling cascade in various non-neuronal systems. The isoform p66SHC does not mediate Ras activation, but is involved in signal transduction pathways that regulate the cellular response to oxidative stress and life span. It acts as a downstream target of the tumor suppressor p53 and is indispensable for the ability of stress-activated p53 to induce elevation of intracellular oxidants, cytochrome c release and apoptosis.

Description

Rabbit polyclonal antibody to SHC

Formulation

Liquid. Neat serum containing 0.05% sodium azide.

Immunogen

Recombinant human SHC portion (aa 1180-1620)

Uniprot ID

P29353

Specificity

Human, Mouse, Hamster

Purification

Serum

Storage

Store at -20°C. Avoid multiple freeze/thaw cycles.

Application

ELISA, IP, WB

Ordering Information

Products	Size	Cat. No.
SHC Polyclonal Antibody	100 µl	A-0494-100

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