

Aryl Hydrocarbon Receptor Polyclonal Antibody

(Catalog No. A-0639)

Description

Rabbit polyclonal antibody to Aryl hydrocarbon receptor

Formulation

Liquid. In TBS containing 0.02% sodium azide.

Immunogen

Recombinant protein corresponding to aa 1-402 of the N-terminal B-1 AhR (Aryl hydrocarbon receptor).

Specificity

Mouse

Purification

Protein A-affinity purified.

Storage

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

Application

ChIP, IP, WB (1:2000)

Ordering Information

Products

Aryl Hydrocarbon Receptor Polyclonal Antibody

Size

25 µg
100 µg

Cat. No.

A-0639-025
A-0639-100

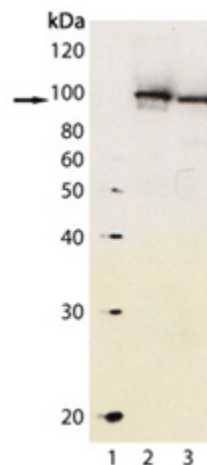


Fig. 1. Western blot analysis of Aryl Hydrocarbon Receptor Polyclonal Antibody: Lane 1: MW marker; Lane 2: Mouse liver extract; Lane 3: Rat liver extract.

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