

96-Well Ras Activation ELISA Kit (Colorimetric)

Our 96-Well Ras Activation ELISA Kit uses the Raf1 RBD (Rho binding domain) bound to a 96-well plate to selectively pull down the active form of Ras from purified or endogenous samples. The captured GTP-Ras is then detected by a pan-Ras antibody and HRP-conjugated secondary antibody.

Ordering Information

[Order Online »](#)

CBI-STA-440	96 assays
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Raf1 RBD detects only the active form of Ras
- Convenient 96-well strip-well format

Handling & Storage

Use/Stability Upon receipt, aliquot and store Raf-1 RBD at -80°C and avoid freeze/thaw. Aliquot and store the anti-pan-Ras Antibody, Secondary Antibody, GTPγS, and GDP components at -20°C and avoid freeze/thaw. Store all other components at 4°C until their expiration dates.

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com