

Human TEAD4/YAP Assay System

This TEAD4/YAP kit product is an all-inclusive assay system that includes in addition to TEAD4/YAP Reporter Cells, two optimized media for use in recovering the cryopreserved cells and for diluting test samples, the reference inhibitor of TEAD Flufenamic acid, Luciferase Detection Reagents, and a cell culture-ready assay plate.

Ordering Information

[Order Online »](#)

| | |
|------------------|-------------|
| NGO-IB16001-0032 | 3x32 assays |
| NGO-IB16001-0096 | 96 wells |
| NGO-IB16002-0384 | 384 wells |

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Transcriptional Enhanced Associate Domain 4 Protein / Yes-Associated Protein 1,
Hippo pathway



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