

# Human TrkA Reporter Assay System

## Ordering Information

[Order Online »](#)

NGO-IB27011-0096	96 wells
NGO-IB27012-0384	384 wells

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Shipping

Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name**

Tropomyosin Receptor Kinase A, NTRK1

**Contents**

This Human Tropomyosin Receptor Kinase A (TrkA; NTRK1) assay kit is an all-inclusive luciferase reporter assay system that includes, in addition to TrkA Reporter Cells, two optimized media for use in recovering the cryopreserved cells and in preparing test samples, reference activator NGF- $\beta$ , Luciferase Detection Reagent, a cell culture-ready assay plate, and a detailed protocol.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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