

NFAT Reporter Assay System

This Nuclear Factor of Activated T cells (NFAT) kit product is an all-inclusive assay system that includes addition to NFAT Reporter Cells, two optimized media for use during cells culture and in diluting test samples, the calcineurin-NFAT pathway activator Ionomycin, Luciferase Detection Reagent, and a cell culture-ready assay plate.

Ordering Information

[Order Online »](#)

NGO-IB18001-0032	3x32 assays
NGO-IB18001-0096	96 wells
NGO-IB18002-0384	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 Nuclear Factor of Activated T Cells



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)