

ROS/SUPEROXIDE DETECTION

Total ROS/Superoxide Detection Kit

ENZ-51010

200 Reactions

HIGHLIGHT

- Directly monitors global levels of reactive oxygen species (ROS), and specifically superoxide, in live cells by fluorescence microscopy or flow cytometry
- Distinguishes between different reactive species, such as hydrogen peroxide, peroxynitrite and hydroxyl radicals
- High sensitivity, specificity and accuracy for live cell studies
- Compatible with major components of tissue culture media (phenol red, FBS and BSA)
- Complete set of reagents, including ROS inducers and scavengers
- Stringently manufactured, to control and eliminate non-specific assay artifacts

Free radicals and other reactive species play seminal roles in many physiological and pathophysiological processes. Once produced within a cell, free radicals can damage a wide variety of cellular constituents, including proteins, lipids and DNA. However, at lower concentrations these very same agents may serve as second messengers in cellular signaling. Information-rich methods are required to quantify the relative levels of various reactive species.

The Total ROS/Superoxide Detection Kit is designed to directly monitor real time reactive oxygen species (ROS) production in live cells using fluorescence microscopy or flow cytometry. The kit includes two fluorescent dyes as major components: Total ROS Detection Reagent (green fluorescent) and specific Superoxide Detection Reagent (orange fluorescent). Through the combination of two specific fluorescent probes, the kit provides a simple and specific assay for the real-time measurement of global levels of reactive oxygen species (ROS), and specifically superoxide in living cells.

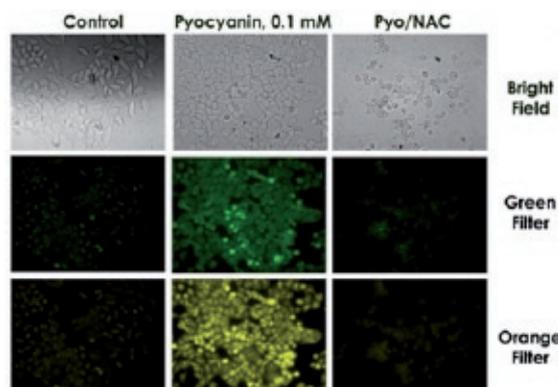


FIGURE 1: Profiling of reactive oxygen species formation by fluorescence microscopy was achieved in HeLa cells loaded with Enzo Life Sciences' Total ROS/Superoxide detection reagents and treated with pyocyanin. General oxidative stress levels were monitored in the green channel, while superoxide production was detected in the orange channel. Pretreatment with NAC, a general ROS inhibitor, prevents formation of ROS.

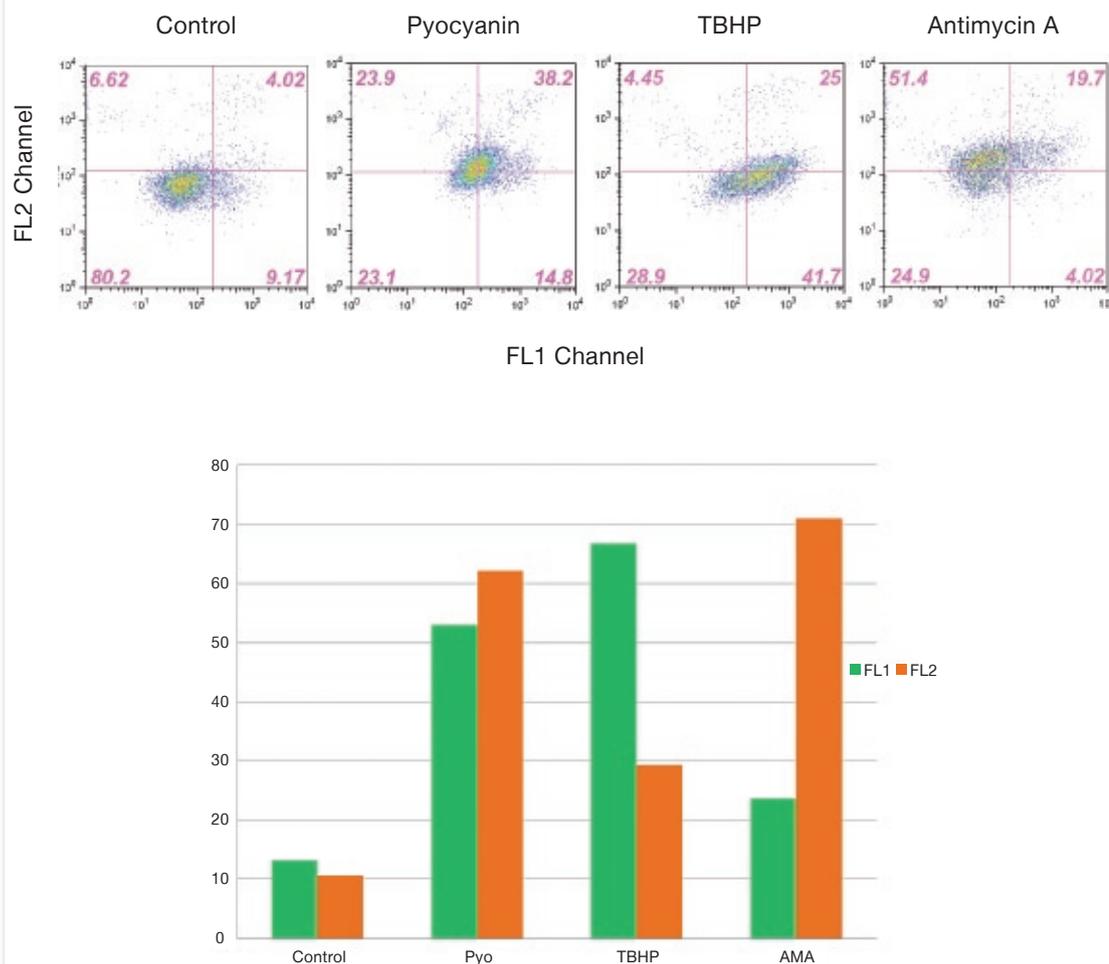


FIGURE 2: Profiling of reactive oxygen species formation by flow cytometry was achieved in Jurkat cells: ROS positive control samples, induced with pyocyanin, yields strong fluorescence signal in both the FL1 and FL2 channels (ROS and superoxide generation); Antimycin A yields strong signal that is detectable in the FL2 channel (superoxide generation); TBHP yields strong signal detectable in the FL1 channel (ROS generation).

Related Products

Product	Prod. No.	Size
ROS/RNS Detection Kit	ENZ-51001	200 Reactions
Red Hydrogen Peroxide Assay Kit	ENZ-51004	500 Reactions
Luminol (Ultra Pure)	ENZ-52354	1 g

Enzo, Alexis and Biomol are trademarks of Enzo Life Sciences, Inc. Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.

Switzerland & Rest of Europe

ENZO LIFE SCIENCES AG

Industriestrasse 17, Postfach
CH-4415 Lausen / Switzerland
Tel. + 41/0 61 926 89 89
Fax + 41/0 61 926 89 79
info-ch@enzolifesciences.com

North/South America

ENZO LIFE SCIENCES INTERNATIONAL, INC.

5120 Butler Pike
Plymouth Meeting, PA 19462-1202
USA
Tel. 1-800-942-0430/(610) 941-0430
Fax (610) 941-9252
info-usa@enzolifesciences.com

Benelux

ENZO LIFE SCIENCES BVBA

Melkerijweg 3
BE-2240 Zandhoven / Belgium
Tel. +32/0 3 466 04 20
Fax +32/0 3 466 04 29
info-be@enzolifesciences.com

Germany

ENZO LIFE SCIENCES GmbH

Marie-Curie-Strasse 8
DE-79539 Lörrach / Germany
Tel. +49/0 7621 5500 526
Toll Free: 0800 6649518
Fax +49/0 7621 5500 527
info-de@enzolifesciences.com

UK & Ireland

ENZO LIFE SCIENCES (UK) LTD.

Palatine House
Matford Court
Exeter EX2 8NL / UK
Tel. 0845 601 1488 (UK customers)
Tel. +44/0 1392 825900 (overseas)
Fax +44/0 1392 825910
info-uk@enzolifesciences.com

For Local Distributors please visit our Website.

incorporating

ALEXIS
BIOCHEMICALS

BIOMOL
INTERNATIONAL

www.enzolifesciences.com