

# PRODUCT INFORMATION

## Oxidative Stress MaxSpec® LC-MS Mixture

Item No. 18701

**Supplied as:** A solution in ethanol (1 µg/ml of each compound)

**Fill Volume:** >1 ml

**Storage:** -20°C

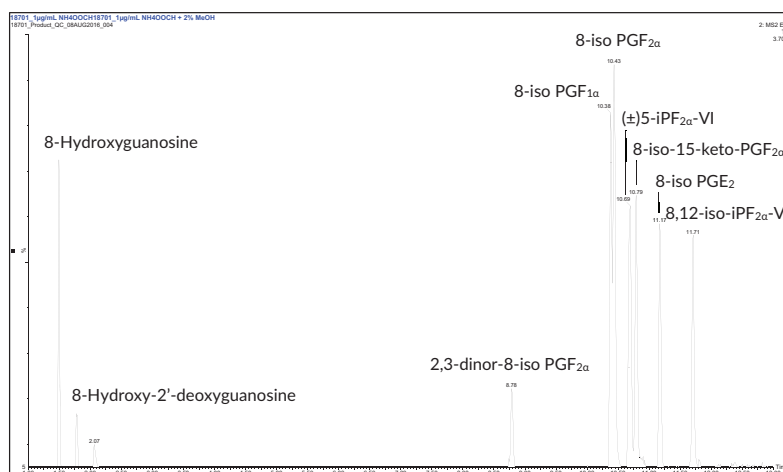
**Stability:** ≥7 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description and Contents

The Oxidative Stress MaxSpec® LC-MS Mixture is a mixture of lipids and nucleic acids produced during oxidative stress in biological systems. The mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation. This product has been designed for direct use in LC-MS applications. The solution may be serially diluted for preparation of calibrators and QC standards and/or used directly as a system suitability standard or tuning standard. After opening, we recommend that the mixture be transferred immediately to a 1 ml glass screw cap vial, to prevent solvent evaporation, and stored at -20°C. The mixture should be discarded after multiple freeze/thaw cycles.

The compounds in this mixture represent the major lipids and nucleic acids that can be detected in serum or urine following oxidative stress. They include the major accepted biomarkers for oxidative stress.



Item Number: 18701		Oxidative Stress MaxSpec® LC-MS Mixture	
Item Number	Item Name	Formula Weight	MS/MS Transition
89300	8-Hydroxyguanosine	299.2	300>168
89320	8-Hydroxy-2'-deoxyguanosine	283.2	284>168
16290	2,3-dinor-8-iso Prostaglandin F <sub>2α</sub>	326.4	325>237
15350	8-iso Prostaglandin F <sub>1α</sub>	356.5	355>293
16350	8-iso Prostaglandin F <sub>2α</sub>	354.5	353>193
16300	(±)5-iPF <sub>2α</sub> -VI	354.5	353>115
16390	8-iso-15-keto Prostaglandin F <sub>2α</sub>	352.5	351>289
14350	8-iso Prostaglandin E <sub>2</sub>	352.5	351>189
16310	8,12-iso-iPF <sub>2α</sub> -VI	354.5	353>115
<b>LC-MS Conditions:</b>			
Mobile Phase A: Water + 0.1% Formic Acid			
Mobile Phase B: Acetonitrile + 0.1% Formic Acid			
Column: Waters BEH C8, 2.1 x 100 mm, 1.7 µm		Flow Rate: 400 µl/min	
0 - 2.0 Minutes: Positive Electrospray Ionization		Full (MS) Scan	
2.0 - 13.0 Minutes: Negative Electrospray Ionization		Full (MS) Scan	

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/17/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM