

# **Oxidative stress markers in high aqueous conditions**

## **Application Note**

Clinical Research

### **Authors**

Agilent Technologies, Inc.

### **Introduction**

Agilent Pursuit PFP columns are ideal for the HPLC analysis of stress markers with water as eluent.



**Agilent Technologies**

## Conditions

Column : Agilent Pursuit PFP, 4.6 x 150 mm, 5  $\mu$ m  
Part No. A3050150X046

Mobile Phase : acetonitrile: 25 mM ammonium acetate pH 3.0 5:95

Flow Rate : 1.0 mL/min

Temperature : Ambient

Detection : 220 nm

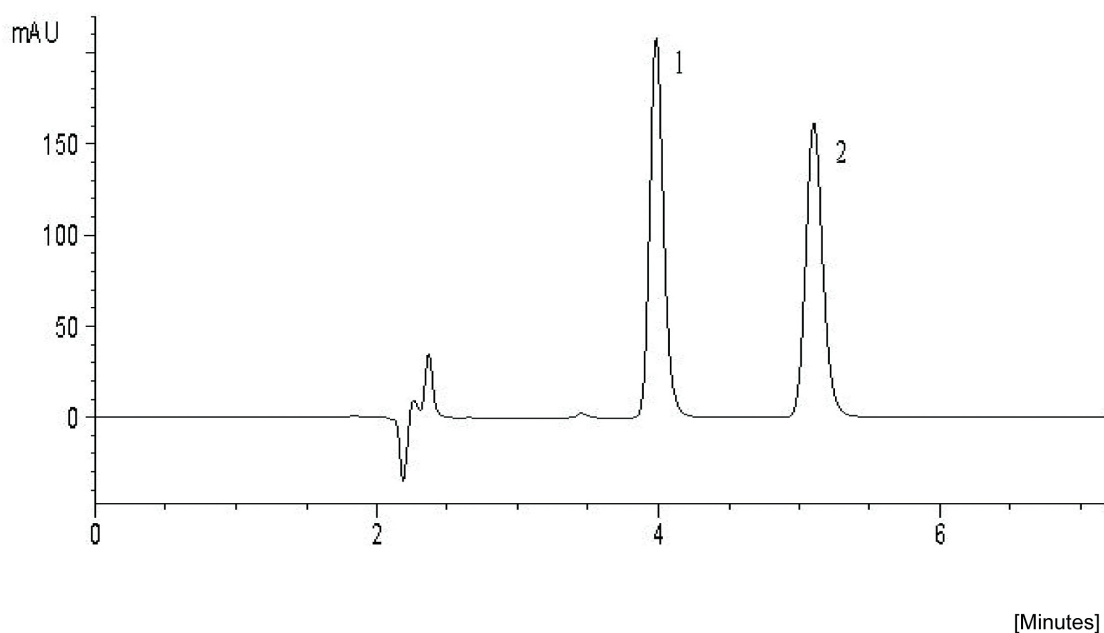
Injection Vol. : 5  $\mu$ L

Sample Concentration : 1 mg/mL all components

Solvent Sample : Water

## Peak identification

1. 2'-deoxyguanosine
2. 8-hydroxy-2'-deoxyguanosine



[www.agilent.com/chem](http://www.agilent.com/chem)

For Research Use Only. Not for use in diagnostic procedures.

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A02451



**Agilent Technologies**