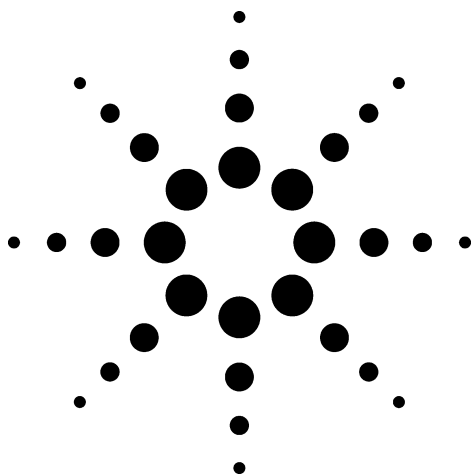




FD&C Colors

Application Food Analysis Robert Ricker

FD&C colorants are commonly used in foods and cosmetics. Propylparaben is a preservative found in the mixture of food colorings used for this analysis. Green food coloring is comprised of yellow # 5 and blue # 1. A combination of ion-pairing and gradient HPLC is used because of the wide chemical differences of compounds. Note the molecular formulas.

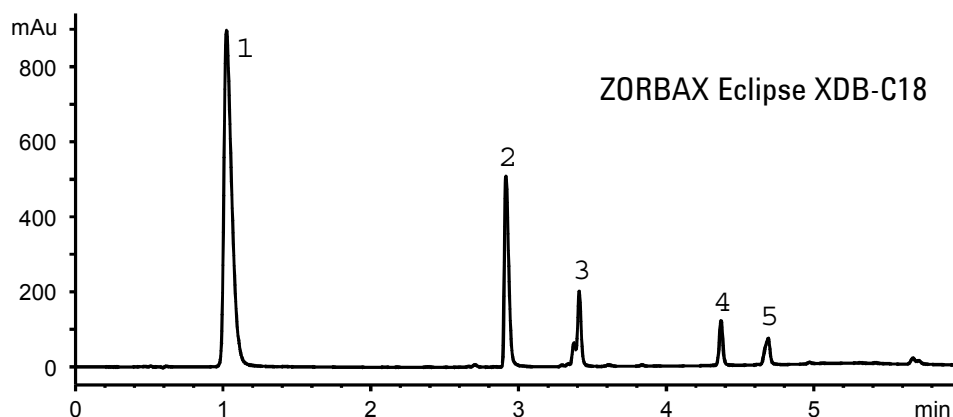
Operating Conditions:

HPLC System: Agilent 1100 with quaternary pump
Column: ZORBAX Eclipse XDB-C18 Rapid-Resolution (3.5 μ m), 4.6 x 50 mm
Agilent Part No. 935967-902
Mobile Phase: A: 0.1% TFA, pH to 4.4 with TEA, B: MeOH
Gradient: 17 to 100% B / 4 min
Detection: UV 254 nm
Flow: 1 mL/ min.
Temperature: ambient

1. Yellow # 5	$C_{16}H_9Na_4Na_3O_9S_2$	MW= 534
2. Red # 40	$C_{18}H_{14}N_2Na_2O_8S_2$	MW= 496
3. Blue # 1	$C_{37}H_{34}N_2Na_2O_9S_3$	MW= 760
4. Propylparaben	$C_{10}H_{12}O_3$	MW= 180
5. Red # 3	$C_{20}H_{14}I_4Na_2O_5$	MW= 878

Highlights

- ZORBAX Eclipse-XDB is designed for extended column life at intermediate pH.
- Reducing column length and particle size simultaneously can:
 - Reduce analysis time
 - Maintain resolution
 - Reduce solvent use



Agilent Technologies

*Robert Ricker is an application chemist
based at Agilent Technologies, Wilmington,
Delaware.*

For more information on our products and
services, visit our website at:
www.agilent.com/chem

Copyright© 2002 Agilent Technologies, Inc.
All Rights Reserved. Reproduction,
adaptation or translation without prior
written permission is prohibited, except as
allowed under the copyright laws.

Agilent shall not be liable for errors
contained herein or for incidental or
consequential damages in connection with
the furnishing, performance, or use of this
material.

Information, descriptions, and specifications
in this publication are subject to change
without notice.

Printed in the USA
April 25, 2002
5988-6355EN



Agilent Technologies