

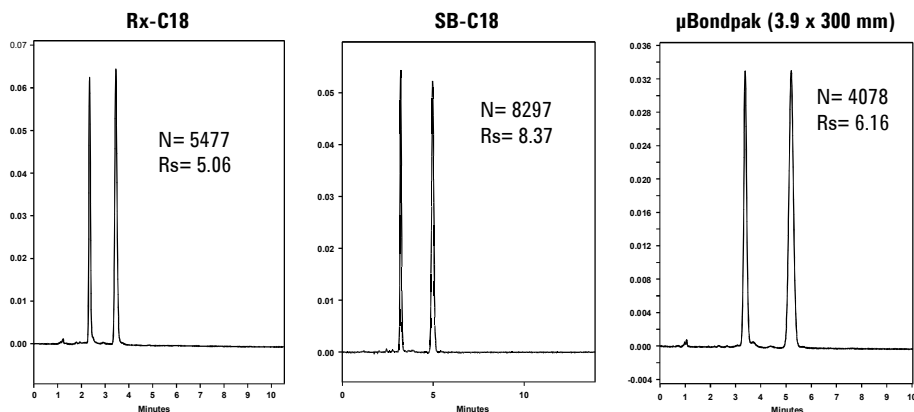
USP Method Triamcinolone and Internal Standard, Hydrocortisone

Application
Pharmaceutical
Robert Ricker

Triamcinolone is a corticosteroid, hence an anti-inflammatory drug, used in combination with other pharmaceuticals in ointments and creams for the treatment of skin eruptions and rashes. It is also used as a single drug product for the treatment of severe bronchial asthma. Triamcinolone was one of the top 30 prescribed generic drugs in 1994.

Highlights

- ZORBAX columns can be substituted for μ Bondapak without compromising the USP criterion that "the resolution factor between triamcinolone and hydrocortisone is not less than 3.0."
- ZORBAX Rx-C18 and SB-C18 both give improved performance demonstrated by higher theoretical plate count as compared to μ Bondapak.
- Highly efficient chromatography and quantitation is easily achieved with both ZORBAX Rx-C18 and SB-C18.



1. Triamcinolone
2. Hydrocortisone

Conditions:
ZORBAX Rx-C18, SB-C18 (4.6 x 250 mm) (Agilent P/N: 880975-902, 880967-902)
Mobile Phase: Water: MeOH, (40:60)
Injection: 10 μ L, 1.5 mL/min, Ambient, Detect. UV(254 nm)



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
5988-6384EN



Agilent Technologies