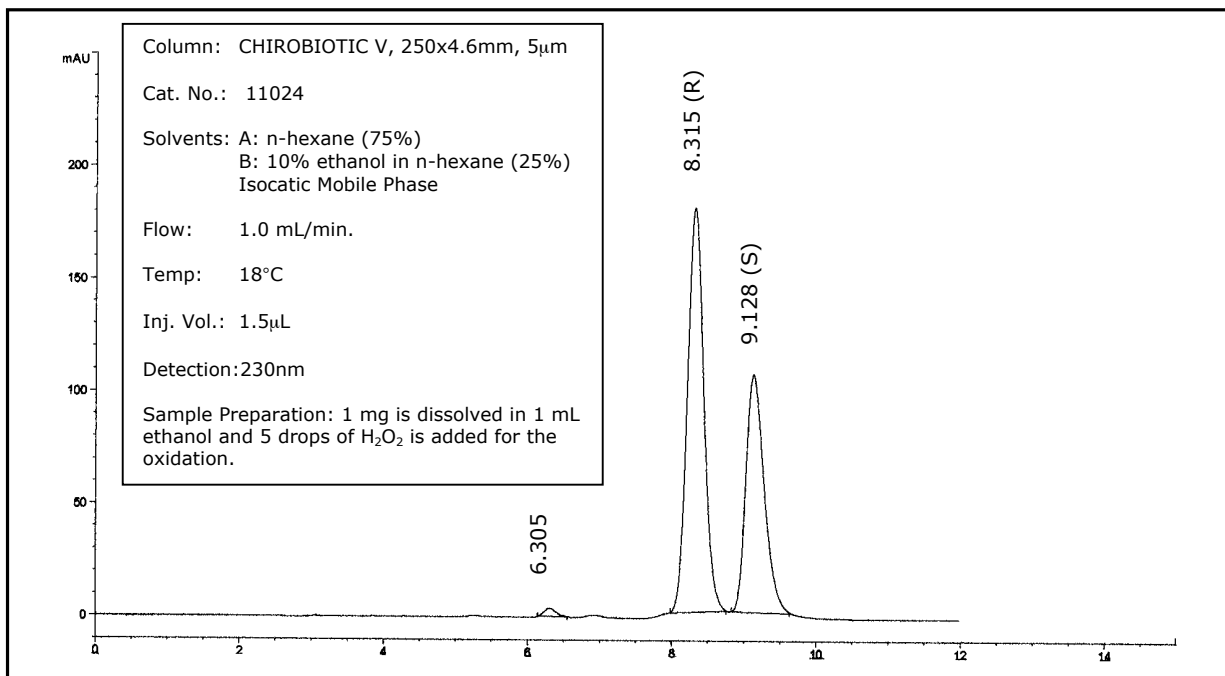
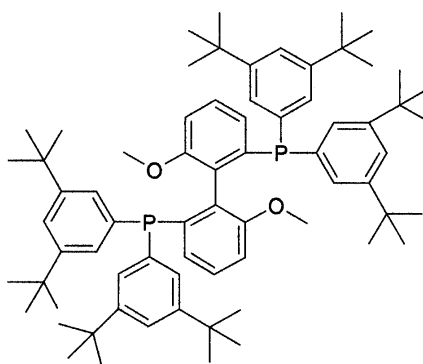


APPLICATION REPORT *astec*

LC001 Phosphine Catalyst

HPLC-Method for the Separation of the Enantiomers of 3,5-tBu-MeOBIPHEP
Using a CHIROBIOTIC V, 250x4.6mm Column

During the synthesis of (6,6'-dimethoxy-2-2'-biphenylene)bis-(bis[3,5-di-tert-butylphenyl]phosphine) (3,5-tBu-MeOBIPHEP), an efficient analytical HPLC chiral separations method was developed using the CHIROBIOTIC V, 250x4.6mm column. Only the bis(oxide) form of the molecule can be separated. Other closely related molecules can be resolved by the same method.



Courtesy of Dr. Marcel Althaus, Hoffman-LaRoche, Basel, Switzerland.

LC001-5/05 ©2005 Advanced Separation Technologies Inc. All rights reserved. Printed in USA.

37 Leslie Court P. O. Box 297 Whippany, NJ 07981 USA
Tel: 973.428.9080 Fax: 973.428.0152 e-mail: info@astecusa.com