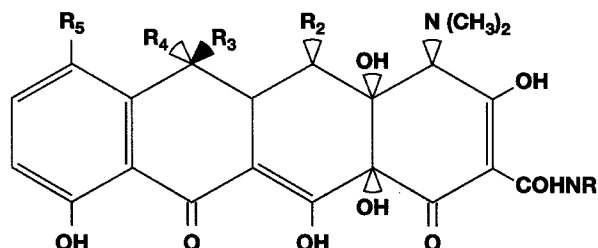


# Separation of Tetracycline Antibiotics by MEKC

Tetracyclines represent a group of antibiotics produced by various *Streptomyces* species. They inhibit protein synthesis in prokaryotes by blocking the binding of aminoacyl tRNAs to ribosomes. Tetracyclines consist of four fused six-membered rings which differ in their substituents (Figure 1). These species can be separated by micellar electrokinetic capillary chromatography (MEKC) using Triton® X-100 as the pseudophase. Because of their large ring structures, tetracyclines can be detected with good sensitivity at 265 nm.



	R <sub>1</sub>	R <sub>2</sub>	R <sub>3</sub>	R <sub>4</sub>	R <sub>5</sub>
<b>Chlortetracycline</b>	H	H	OH	CH <sub>3</sub>	Cl
<b>Oxytetracycline</b>	H	OH	OH	CH <sub>3</sub>	H
<b>Tetracycline</b>	H	H	OH	CH <sub>3</sub>	H
<b>Epitetracycline</b>	H		CH <sub>3</sub>	OH	

## Analysis Conditions

Instrument	BioFocus® 3000 TC system
Capillary	24 cm x 25 µm, coated with linear polyacrylamide
Run buffer	200 mM sodium phosphate + 0.12% reduced Triton X-100 (pH 2.2)
Purge protocol	60 seconds, with run buffer
Injection	pressure at 10 psi * second
Polarity	positive to negative
Voltage	10 kV (observed current 36 µA)
Detection	265 nm
Cartridge temperature	20 °C
Autosampler temperature	20 °C

Fig. 1. Structures of tetracycline antibiotics.

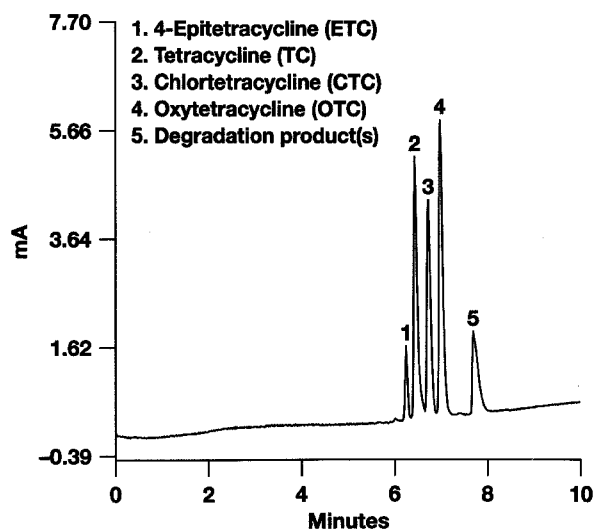


Fig. 2. Separation of tetracyclines by MEKC.

Triton is a trademark of Rohm and Haas.

**Bio-Rad  
Laboratories**

ISO 9001 registered

www.bio-rad.com

Life Science  
Group

U.S. (800) 4BIORAD Australia 02-9914-2800 Austria (1)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771 China (86-10) 2046622  
Denmark 39 17 9947 Finland 90 804 2200 France (1) 43 90 46 90 Germany 089 31884-0 Hong Kong 7893300 India 91-11-461-0103  
Israel 03 951 4127 Italy 02-21609.1 Japan 03-5811-6270 The Netherlands 0313 18-540666 New Zealand 09-443 3099  
Singapore (65) 272-9877 Spain (91) 661 70 85 Sweden 46 (0) 8 627 50 00 Switzerland 01-809 55 55 United Kingdom 0800 181134