

**EUROPE**

LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM • Tel.: +32 4 372 74 00  
• Fax: +32 4 372 75 00 Toll-free: + 800 666 00 123 • info@eurogentec.com •  
[www.eurogentec.com](http://www.eurogentec.com)

**NORTH AMERICA**

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA • Tel.: +1-510-791-9560  
Toll-free: +1 800-452-5530 • Fax: +1 510-791-9572 • [service@anaspec.com](mailto:service@anaspec.com)  
[www.anaspec.com](http://www.anaspec.com)



Experience true partnership

Eurogentec products are sold for research or laboratory use only and are not to be administered to humans or used for medical diagnostics

## Technical Data Sheet

# BioMarker™ CST 524/420

MW-0200-SP

[250µL]

### Description:

Biomarkers are optimal for reading DNA fragments ranging from 50 to 3000 bp.

BioMarker™ CST 524/420 contains linear double stranded DNA bands of 1200, 800, 524, 420, 373, 200, 100, and 50 base pairs.

The concentration of each band is at 50ng/µL of BioMarker™ CST 524/420 applied.

### Shipping conditions:

The BioMarker™ CST 524/420 is shipped at room temperature

### Storage conditions:

The BioMarker™ CST 524/420 must be stored at 4°C

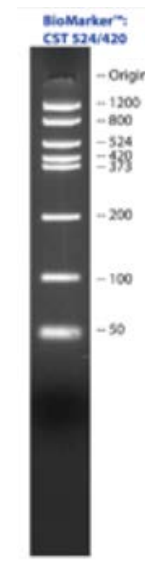


Figure 1: BioMarker™ CST 524/420 Profil on a 4% NuSieve © 3:1 Plus Agarose gels

**For further information please contact our Customer Help Desk:**

**For Europe:**

E-mail: [info@eurogentec.com](mailto:info@eurogentec.com)  
Tel.: +32 4 372 76 65 • Toll-free: + 800 666 00 123

**For USA:**

E-mail: [service@anaspec.com](mailto:service@anaspec.com)  
Tel.: +1-510-791-9560 • Toll-free: +1-800-452-5530