

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

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
• www.anaspec.com

Technical Data Sheet

iD Graphite Electrode **ID-STETD1-001 • ID-BLETD1-001**

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

Product

Cat #	Description	Size
ID-STETD1-001	iD Stain Graphite Electrode	1 unit
ID-BLETD1-001	iD Blot Graphite Electrode	1 unit

Description

The iD Stain or Blot Graphite Electrode made of powered carbon and clay, serves as the anode replacement of the iD Stain Gel Device or iD Blot Device respectively.

Treated with a special process, each unit of the Graphite Electrode possesses excellent electrical conduction, stable electrochemical structure, adequate mechanical strength, and good exhaust function.

Specifications

Dimensions: 100 mm (l) × 92 mm (w) × 8.8 mm (t)

Weight: 206 g

Materials: Powdered carbon, Clay, Stainless steel

For further information please contact our Customer Help Desk:

For Europe:

E-mail: info@eurogentec.com

Tel: +32 4 372 76 65

For USA:

E-mail: service@anaspec.com

Tel: +1 510 791 9560