

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures.

CometSlide™ Rack System

Catalog #: 4252-040-02

Description:

The CometSlide™ Rack System was designed to accommodate standard CometSlides™ and FLARE™ slides as wells as CometSlide™ HT for use in CometAssay™ and FLARE™ assays.

Contents:

Each CometSlide™ Rack System is composed of a durable plastic rack that holds 20 slides, and a translucent plastic reservoir with lid.

Instructions:

For CometSlide™ (1" x 3"), add 200 ml of buffer to reservoir; rack slides; submerge slides in buffer; and incubate as directed.

For CometSlide HT (2" x 3"), add 350 ml of buffer to reservoir; rack slides; submerge slides in buffer; and incubate as directed.

Product Care:

A mild laboratory detergent may be used, if desired, followed by thorough rinsing with deionized water. The system may be sprayed down with 70% Ethanol for decontamination. Rinse thoroughly with deionized water before each use. **Do not Autoclave! Exposure to ultraviolet light for extended periods of time may compromise integrity of product. Not compatible with acetone or xylenes.**

Storage:

Store at Room Temperature.

©2005, Trevigen, Inc. CometSlide and CometAssay are trademarks and Trevigen is a registered trademark of Trevigen, Inc.

TREVIGEN, INC.

8405 Helgerman Court, Gaithersburg, MD 20877 USA
Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800
Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures.

CometSlide™ Rack System

Catalog #: 4252-040-02

Description:

The CometSlide™ Rack System was designed to accommodate standard CometSlides™ and FLARE™ slides as wells as CometSlide™ HT for use in CometAssay™ and FLARE™ assays.

Contents:

Each CometSlide™ Rack System is composed of a durable plastic rack that holds 20 slides, and a translucent plastic reservoir with lid.

Instructions:

For CometSlide™ (1" x 3"), add 200 ml of buffer to reservoir; rack slides; submerge slides in buffer; and incubate as directed.

For CometSlide HT (2" x 3"), add 350 ml of buffer to reservoir; rack slides; submerge slides in buffer; and incubate as directed.

Product Care:

A mild laboratory detergent may be used, if desired, followed by thorough rinsing with deionized water. The system may be sprayed down with 70% Ethanol for decontamination. Rinse thoroughly with deionized water before each use. **Do not Autoclave! Exposure to ultraviolet light for extended periods of time may compromise integrity of product. Not compatible with acetone or xylenes.**

Storage:

Store at Room Temperature.

©2005, Trevigen, Inc. CometSlide and CometAssay are trademarks and Trevigen is a registered trademark of Trevigen, Inc.

TREVIGEN, INC.

8405 Helgerman Court, Gaithersburg, MD 20877 USA
Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800
Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com

CometSlide™ Rack System
Catalog #: 4252-040-02
Storage: Room Temperature
TREVIGEN®
1-800-873-8443

CometSlide™ Rack System
Catalog #: 4252-040-02
Storage: Room Temperature
TREVIGEN®
1-800-873-8443