

# Getting Started with the Ultraviolet Sample Tray

An ultraviolet sample tray (catalog #170-8271) is available for the Gel Doc™ EZ imaging system. To get started using the ultraviolet tray, perform the following steps.

## Clean the Tray

Clean the tray with standard laboratory detergents or mild solvents such as ethanol or methanol. Use lint free tissue to wipe it dry. Dust particles or lint on a sample tray can glow under UV illumination.

Caution: Do not clean the ultraviolet tray in a dishwasher. Doing so will damage the tray.

## Turn on the System

Turn on the imager and start Image Lab™ software.

## Insert the Tray in the Imager

Insert the tray into the Gel Doc EZ imager and close the door. A wizard appears on the computer to guide you through the steps of initializing the tray.

For more information about initializing trays, refer to the Gel Doc™ EZ System Installation Guide.

## Supported Applications

The ultraviolet sample tray supports the following applications:

- Ethidium bromide
- SYBR® Green
- SYBR® Safe
- SYBR® Gold
- GelGreen
- GelRed
- Flamingo™ fluorescent gel stain
- Oriole™ fluorescent gel stain
- Coomassie Fluor Orange
- SYPRO Ruby
- Krypton

## Legal Notices

Coomassie is a trademark of BASF Aktiengesellschaft. GelRed and GelGreen are trademarks of Biotium, Inc. Krypton is a trademark of Thermo Fisher Scientific Inc. SYBR® and SYPRO are trademarks of Invitrogen Corporation.



## 紫外线样本托盘简介

用于 Gel Doc™ EZ 成像系统的紫外线样本托盘可获得（目录编号：170-8273）。  
要开始使用蓝色样本托盘，请执行以下步骤。

### 清洁托盘

使用标准的实验室清洁剂或温和的溶剂（如乙醇或甲醇）清洁托盘。使用无尘软布将其擦干。样本托盘上的微尘或绒毛会在紫外线照射下发光。

注意事项：请勿在洗碗机中清洁紫外线托盘。这样做会损坏托盘。

### 打开系统

打开成像仪并启动 Image Lab™ 软件。

### 将托盘插入成像仪中

将托盘插入 Gel Doc EZ 成像仪并关上托盘门。

计算机上将出现一个向导，引导您完成初始化托盘的步骤。

有关初始化托盘的更多信息，请参阅《Gel Doc™ EZ 系统安装指南》。

### 支持的应用

紫外线样本托盘支持以下应用：

- Ethidium bromide
- SYBR® Green
- SYBR® Safe
- SYBR® Gold
- GelGreen
- GelRed
- Flamingo™
- Oriole™
- Coomassie Fluor Orange
- SYPRO Ruby
- Krypton

### 法律声明

Coomassie 是 BASF Aktiengesellschaft 的商标。GelRed 和 GelGreen 是 Biotium, Inc 的商标。Krypton 是 Thermo Fisher Scientific Inc 的商标。SYBR® 和 SYPRO Invitrogen Corporation 的商标。