

# Gel Alignment Template Kit

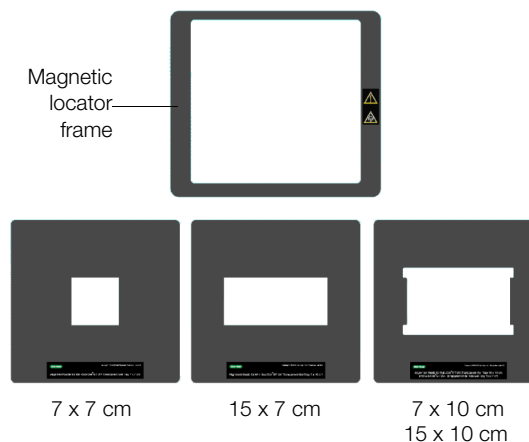
The Bio-Rad gel alignment template kit allows four sizes of standard agarose gels to be centered quickly and easily, as well as ensuring consistent placement of gels over time. The gel alignment template kit is offered for the convenience of Gel Doc™ XR and ChemiDoc™ XRS camera system users.

## Gel Alignment Template Kit Contents

The gel alignment template kit (catalog #170-8184) contains:

- Magnetic locator frame
- Alignment guide for Sub-Cell® GT UV-transparent mini-gel tray, 7 x 7 cm
- Alignment guide for Sub-Cell GT UV-transparent wide mini-gel tray, 15 x 7 cm
- Alignment guide for Sub-Cell GT UV-transparent mini-gel tray, 7 x 10 cm and Sub-Cell GT UV-transparent gel tray, 15 x 10 cm

The gel alignment templates can be used with the XcitaBlue™ UV to blue conversion screen (catalog #170-8183). The templates fit exactly into the XcitaBlue conversion screen frame.



## Gel Alignment Template Kit Instructions

1. Place the locator frame over the transilluminator, with the magnetic side down. Match the corners of the magnetic locator frame with the edges of the transilluminator image area. The UV symbol on the frame will be in the same orientation as the UV symbol on the imager.
2. Place the gel alignment template that matches the size of your sample tray or agarose gel into the magnetic locator frame.
3. Place your gel or gel tray into the open area of the template.

