

PlateLoc Sealer Customer Notice

Potential adjustment required before using Peelable Aluminum RT

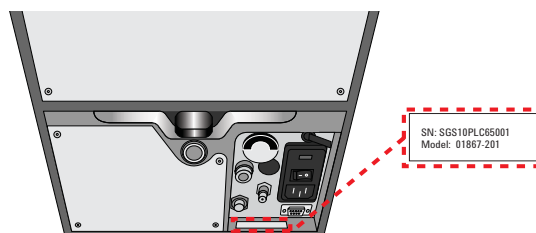
Before using the Peelable Aluminum RT heat seal or sample roll (part numbers 24214-001 and 24214-081, respectively), determine if your PlateLoc Sealer requires an adjustment to ensure compatibility.

CAUTION The Peelable Aluminum RT is thicker than previous Agilent Technologies heat seal material and can become wedged within the material-feeding channel of older PlateLoc Sealers. The material could be difficult to remove and the device could require service. Before using the thicker material, you should determine if your device requires an adjustment.

Determining if your PlateLoc Sealer requires an adjustment

PlateLoc Sealers manufactured before week 51 in December 2010 might require an adjustment to use the thicker Peelable Aluminum RT seal material. To determine if your device should be evaluated for compatibility, check the device serial number. You can locate the serial number (SN) label at the back of the device below the connection panel.

Figure SN label on PlateLoc Sealer (back view)



Does the serial number (SN) start with the following prefix?	SN description or example	Evaluation required
SGS10	The third and fourth digits from the end are 51 or 52, for example: SGS10PLC65101 SGS10PLC45201	No
	The third and fourth digits from the end are 50 or lower, for example: SGS10PLC65001 , SGS10PLC64901 , and so forth	Yes See Adjustment details below.
Any prefix other than SGS	For example, USS2009180002	Yes See Adjustment details below.

PlateLoc Sealer Customer Notice

A serial number that starts with an SGS11 prefix or greater indicates that the PlateLoc Sealer was manufactured later than 2010. Such devices have already been adjusted to ensure compatibility with the Peelable Aluminum RT.

Adjustment details

If your PlateLoc Sealer requires evaluation for compatibility with the Peelable Aluminum RT, obtain a sample roll of the heat seal and schedule the evaluation during an Agilent Technologies instrument service visit (for example, annual Preventive Maintenance). During a scheduled service visit, the Agilent Field Service Engineer can make the necessary PlateLoc Sealer adjustment without an additional charge. Typical adjustments include the Seal Entry Slot and the Cutter Blade. After the adjustments are made, the instrument will not require further adjustment when alternating between the Agilent heat seal material of different thicknesses.

Contact information

For more information, contact one of the following:

Technical Support: 1.800.979.4811 or +1.408.345.8011
service.automation@agilent.com

Customer Service: 1.866.428.9811 or +1.408.345.8356
orders.automation@agilent.com

European Service: +44 (0)1763853638
euroservice.automation@agilent.com

Web: www.agilent.com/lifesciences/automation