

EVOWare Manual Instructions

Dear Freedom EVO user,

To ensure your instrument continues to perform optimally, we recommend that you update the **Tip Offset** value in your EVOware software (see Figure 1) using the new settings shown in Table 2. This update applies to the part numbers listed in Table 1.

Entering these new values correctly is important to avoid workflow disruptions.

If you have any questions or need help with the update process, please do not hesitate to [contact us](#).

Thank you for your attention to this important update.

Table 1. Part numbers.

Tip size	Refill tray format	Applicable part numbers
1000 µl	SLAS	30201342, 30199162
	Hanging	10612554, 10612555, 10612513, 30064860, 30064861
1000 µl WB	SLAS	30201344
	Hanging	30201343
200 µl	SLAS	30201340, 30201341
	Hanging	10612552, 10612553, 30064848, 30064849
50 µl	SLAS	30201338, 30201339
	Hanging	30200713, 30200712

Table 2. Tip Offset values.

Tip size	New Tip Offset values
1000µl	332
200µl	-43
50µl	-50

Figure 1. Labware Editor.

