

Dharmacon™ CRISPR RNA Configurator User Guide has been replaced. Please see the new version of the [CRISPR Design Tool - User Guide](#).