



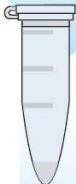
## MIX

Mix RNA with kit contents

8-50  $\mu$ l RNA

0.1x volume Reaction Buffer

1  $\mu$ l HL-dsDNase



## HEAT

Incubate

10 minutes, 37°C

5 minutes, 55°C



## RUN

Run RT of your choice  
in the same tube