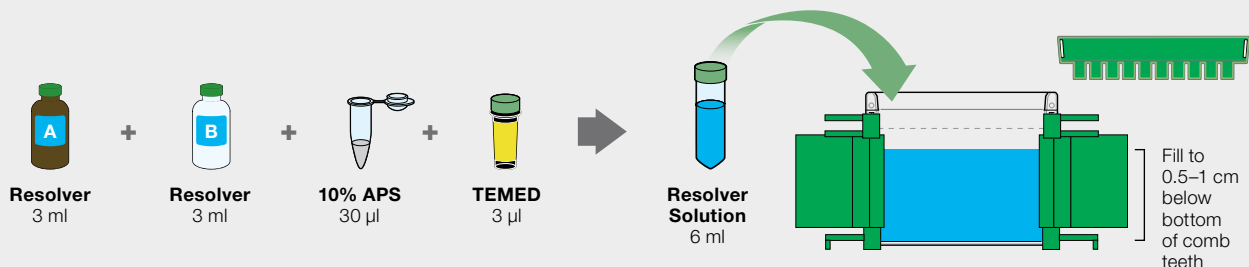


# How to Cast Your Gels Fast

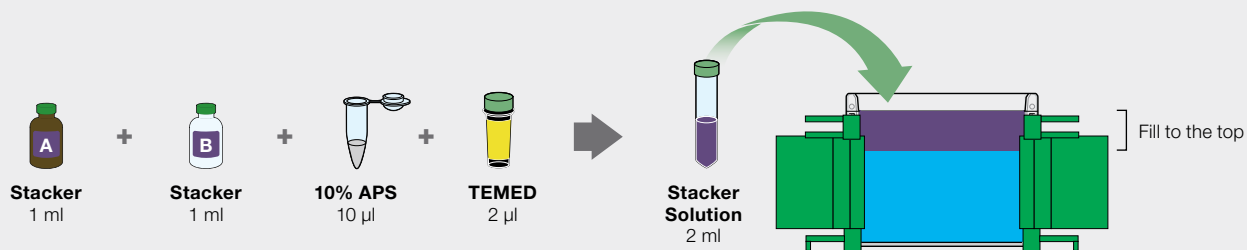
This quick guide illustrates how to cast an individual gel with Bio-Rad's TGX™ and TGX Stain-Free™ FastCast™ acrylamide kits using 1.0 mm Bio-Rad glass plates.

To cast multiple gels using these kits with 1.0 mm Bio-Rad glass plates, multiply the amounts below by the number of gels desired. For information on using other casting setups, please see Table 2 in the instruction manual.

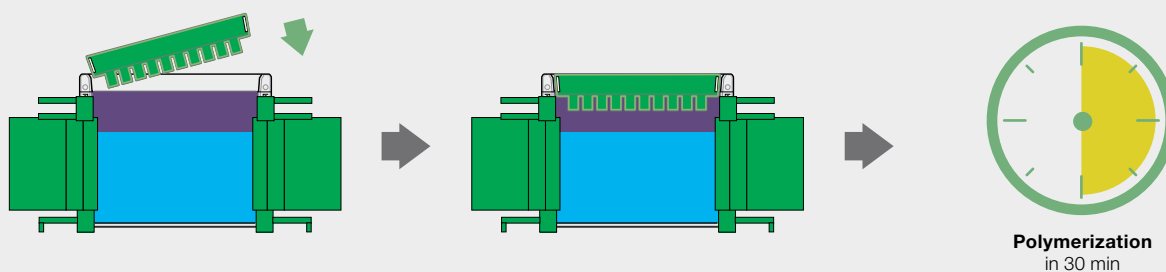
## 1 Prepare Resolver Solution



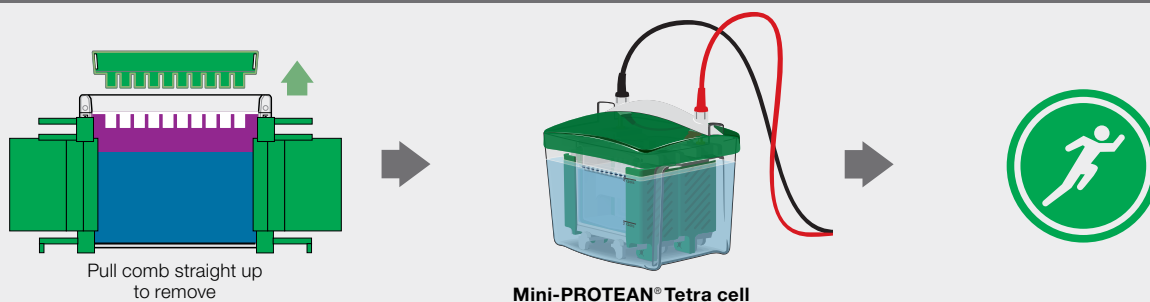
## 2 Prepare Stacker Solution



## 3 Insert Comb and Allow Polymerization



## 4 Run Gel



**BIO-RAD**