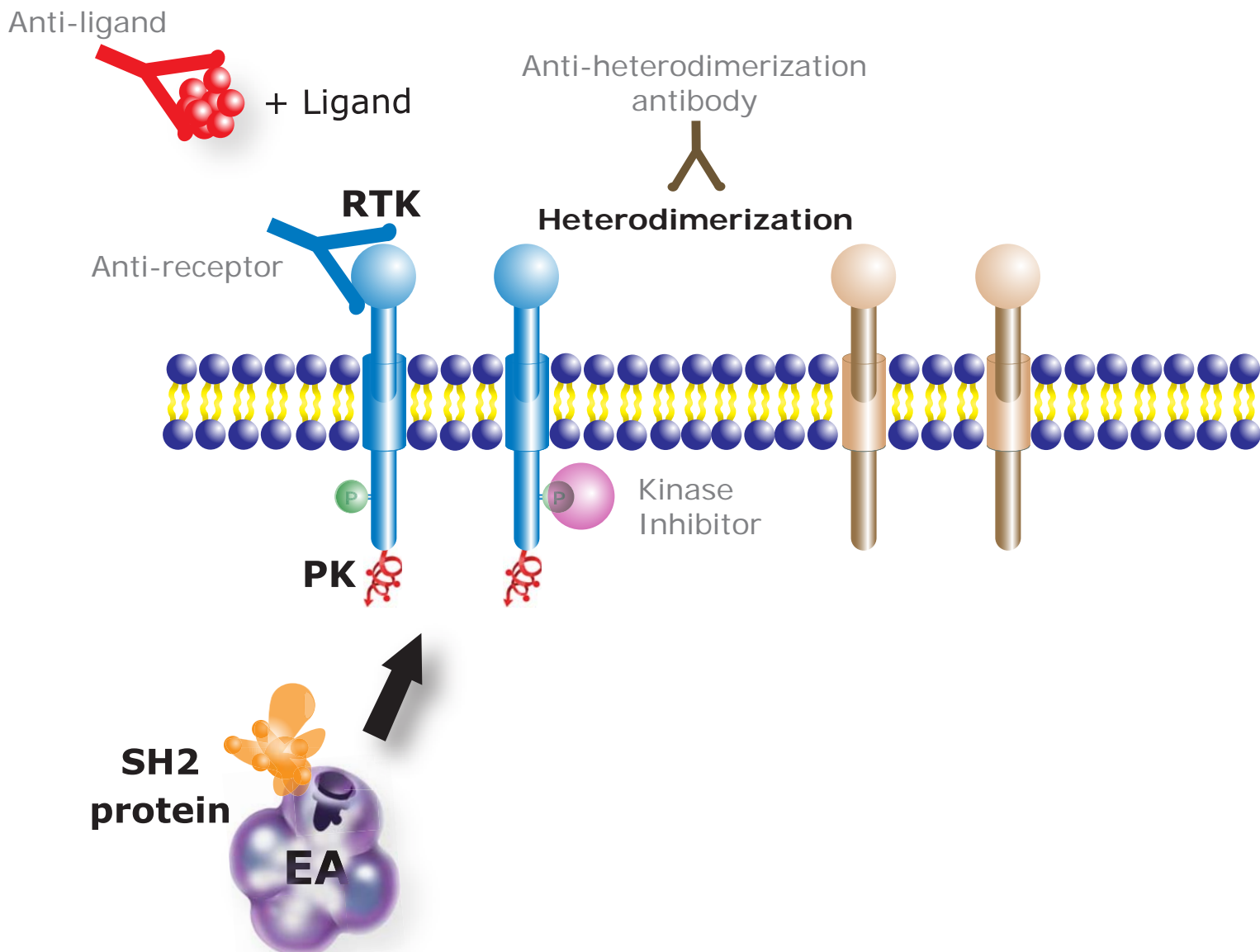


Novel Platform for therapeutic Antibody testing



Use of a full length receptor protein allows identification of novel antibodies in PathHunter assays.