

ThINQ!™ Activity



Help your students design and plan inquiry investigations

Use this design and planning worksheet to help your students engage in inquiry investigations.

Each section guides students through the process of designing and planning an experiment just like a real scientist. This can be used as a homework assignment or an in-class exercise. Review these worksheets to assess the strengths and weaknesses of your students' understanding of content and the scientific process. It will also help you plan and prepare materials and equipment needed by your students to complete their inquiry investigations. This worksheet is a perfect complement to Bio-Rad's ThINQ! Investigation Kits. Visit bio-rad.com/thinq for more information.



Title:

Authors:

Hypothesis (1–2 sentences):

Background (3–4 sentences):

Questions:

Variables (dependent and independent):

Constants:

Safety Notes:

Proposed Procedure:

Written steps	Drawn steps

What data will be collected?

How will the data be analyzed? Sketch out what the data will look like if they support the hypothesis.