

Name:
Grade Level/Subject Taught:
Date:

CureCancer Reflection

- 1. Was this online unit accessible and helpful?*
- 2. How could you use this unit in your classroom?*
- 3. Do you think this unit would help your ELL students understand and participate in Bio-Rad's Secrets of the Rain Forest kit? Why?*
- 4. What improvements would you make? Why?*
- 5. Anything else?*