

- Wash away unbound dye-labeled lysate proteins and quantify captured proteins

