

Building a PathHunter® Protein translocation cell line

