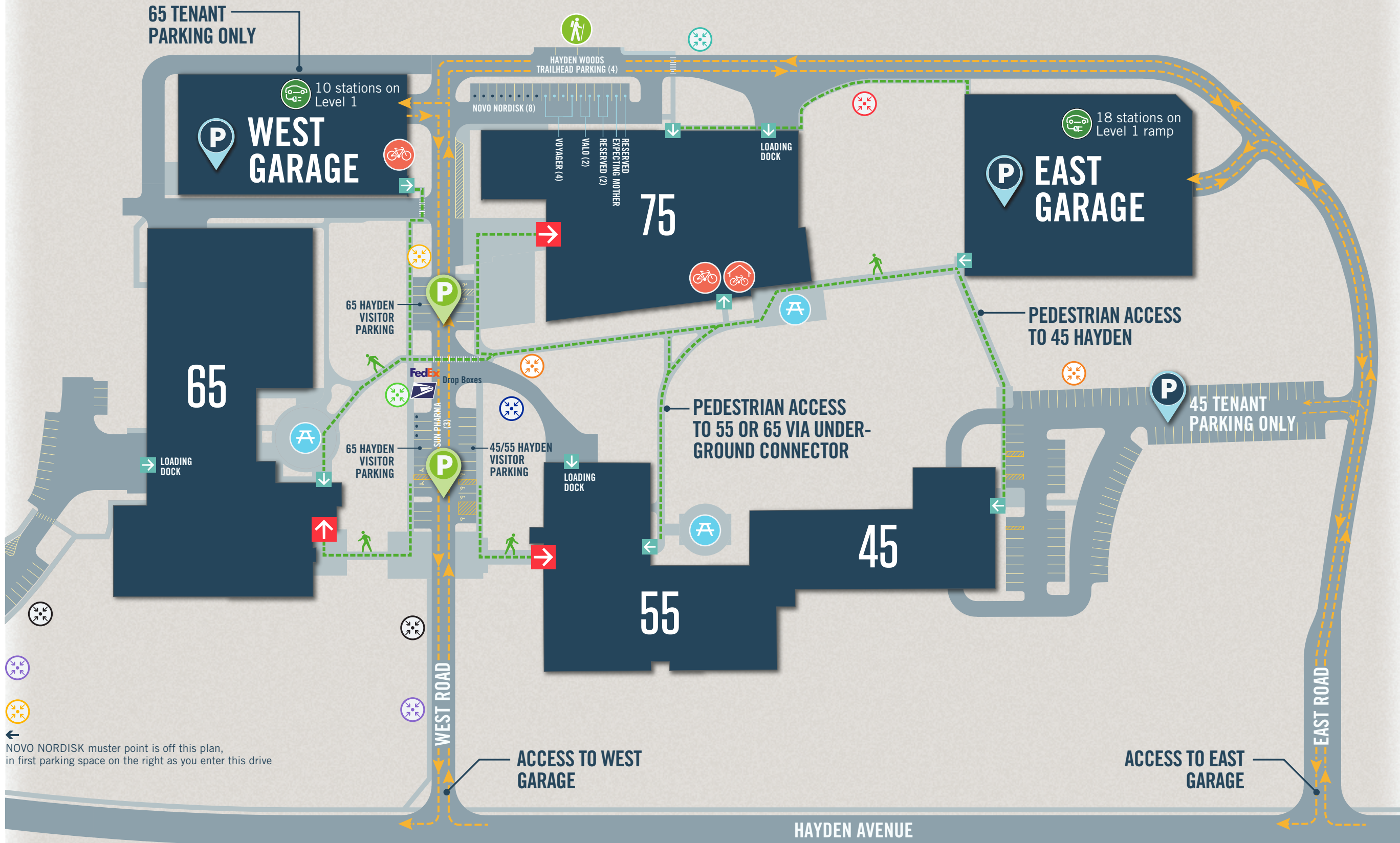


PARKING ACCESS PLAN

HAYDEN RESEARCH CAMPUS



NOVO NORDISK muster point is off this plan, in first parking space on the right as you enter this drive

MUSTER POINT LOCATIONS

- VOYAGER
- NOVO NORDISK 65 HAYDEN
- NOVO NORDISK 75 HAYDEN
- SUN PHARMA
- ALEXION
- ACCENT
- VALO
- TAKEDA

MAP KEY

VISITOR / TENANT

- TENANT PARKING
- VISITOR PARKING
- PEDESTRIAN CIRCULATION
- TENANT VEHICULAR CIRCULATION
- OUTDOOR SEATING
- MAIN ENTRANCE
- ENTRY/EXIT
- ELECTRIC VEHICLE CHARGING
- TRAILHEAD PARKING
- BIKE STORAGE ROOM
- BIKE RACKS
- MUSTER POINT

KSP King Street Properties
Placemaking. For Science.