

# GIPZ<sup>TM</sup> Lentiviral shRNA— glycerol stocks

Applies to all GIPZ Lentiviral shRNA products provided in tubes and microtiter plates as glycerol stocks.

## Product description

Dharmacon<sup>TM</sup> GIPZ<sup>TM</sup> Lentiviral shRNA plasmid vectors are supplied in transformed bacterial cultures grown in the presence of glycerol for storage, either in tubes or microtiter plates.

## Quality control

All cultures are checked for growth before shipment.

## Conclusion

Individual constructs are shipped as bacterial cultures of *E. coli* (PrimePlus) in 2x LB broth (low salt) medium with 8% glycerol, 100 µg/mL carbenicillin and 25 µg/mL zeocin. Individual constructs are shipped on wet ice. Collections are shipped in 96-well plate format on dry ice. Individual constructs and collections should be stored at –80 °C. **To allow any CO<sub>2</sub> that may have dissolved into the media from the dry ice in shipping to dissipate, please store constructs at –80 °C for at least 48 hours before thawing.**

## Important safety note

Please follow the safety guidelines for use and production of vector-based lentivirus as set by your institution's biosafety committee.

- For glycerol stocks of *E. coli* containing lentiviral plasmids, BSL1 guidelines should be followed
- For handling and use of lentiviral products to produce lentiviral particles, BSL2 or BSL2+ guidelines should be followed
- For handling and use of lentiviral particle products, BSL2 or BSL2+ guidelines should be followed

**Please note that GIPZ vectors are not compatible with third generation packaging systems such as ViraPower<sup>TM</sup>. We recommend the Trans-Lentiviral shRNA packaging system for use with our vectors.**

## If you have any questions contact

**t** +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

**f** +44 (0)1223 655 581

**w** [horizondiscovery.com/contact-us](http://horizondiscovery.com/contact-us) **or** [dharmacon.horizondiscovery.com/service-and-support](http://dharmacon.horizondiscovery.com/service-and-support)

**Horizon Discovery**, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom