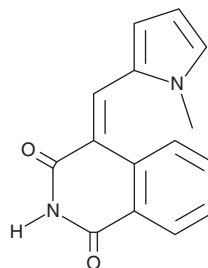


# PRODUCT INFORMATION



**BYK204165**  
Item No. 21430

**CAS Registry No.:** 1104546-89-5  
**Formal Name:** 4-[(1-methyl-1H-pyrrol-2-yl)methylene]-  
1,3(2H,4H)-isoquinolinedione  
**Synonym:** PARP Inhibitor XIV  
**MF:** C<sub>15</sub>H<sub>12</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 252.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 427 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

BYK204165 is supplied as a crystalline solid. A stock solution may be made by dissolving the BYK204165 in the solvent of choice, which should be purged with an inert gas. BYK204165 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BYK204165 in these solvents is approximately 10 mg/ml.

## Description

BYK204165 is a selective inhibitor of poly(ADP-ribose) polymerase 1 (PARP1; IC<sub>50</sub> = 44.67 nM for human recombinant PARP1 in an enzyme assay).<sup>1</sup> It is selective for PARP1 over PARP2 (IC<sub>50</sub> = 4,168 nM for murine PARP2 in an enzyme assay). BYK204165 inhibits hydrogen peroxide-activated PARP in A549 human lung epithelial, C4I human cervical, and H9c2 rat cardiac myoblast cells with IC<sub>50</sub> values of 229.09, 1,778.28, and 123.03 nM, respectively.

## Reference

1. Eltze, T., Boer, R., Wagner, T., *et al.* Imidazoquinolinone, imidazopyridine, and isoquinolindione derivatives as novel and potent inhibitors of the poly(ADP-ribose) polymerase (PARP): A comparison with standard PARP inhibitors. *Mol Pharmacol.* **74**(6), 1587-1598 (2008).

### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/23/2022

## CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)